

Protein Thermal Shift™ studies

USER GUIDE

Using Protein Thermal Shift™ reagents and Protein Thermal Shift™
Software v1.4

Catalog Numbers 4461146, 4462263, 4466037, 4466038

Publication Number MAN0025600

Revision B.0



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN0025600

Revision	Date	Description
B.0	4 October 2022	<ul style="list-style-type: none">Added important note with instructions to not set the temperature of the instrument below room temperature when you load and start a run.Added instructions to prepare files with QuantStudio™ Design and Analysis Software v2, QuantStudio™ Design and Analysis Desktop Software, and QuantStudio™ Real-Time PCR Software.Corrected the recommended filter settings.
A.0	12 July 2021	<p>Pub No. MAN0025600 Rev. A.0 replaces Pub No. 4461808 Rev. A, and includes the following changes from the previous version:</p> <ul style="list-style-type: none">Quantity sufficient for reactions is now 1000

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

NOTICE TO PURCHASER: DISCLAIMER OF LICENSE: Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Thermo Fisher Scientific, either expressly, or by estoppel.

TRADEMARKS: Microsoft and Windows are trademarks of Microsoft Corp. Intel and Pentium are trademarks of Intel Corporation. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2022 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■ CHAPTER 1 About this guide	9
Purpose	9
User attention words	9
■ CHAPTER 2 Getting started with a Protein Thermal Shift™ study	10
Product information	11
Purpose of the starter kit	11
Starter kit contents and storage	11
Purpose of the dye kit	11
Dye kit contents and storage	11
Materials required but not included	12
General guidelines	12
Protein considerations	12
Experimental conditions	13
Plate layout	13
Replicates	13
Controls	13
Supported real-time PCR systems	14
■ CHAPTER 3 Create and set up an experiment file for the instrument run	15
Set up a plate file for high resolution melt analysis in QuantStudio™ Design and Analysis Software v2 (v2.x)	15
Set up a melt curve run in QuantStudio™ Design and Analysis Desktop Software (v1.x) .	18
Design an experiment in QuantStudio™ Real-Time PCR Software (v1.x)	22
Design an experiment in QuantStudio™ 12K Flex Software	25
Create and set up an experiment file for the ViiA™ 7 Real-Time PCR System	28
Create and set up an experiment file for the StepOne™ Real-Time PCR System or StepOnePlus™ Real-Time PCR System	31
Create and set up an experiment file for the 7500 Real-Time PCR System or 7500 Fast Real-Time PCR System	34

■ CHAPTER 4 Prepare and run the reactions, then review results	36
About the example study protein melt reactions	36
Required materials	36
Prepare the protein melt reactions	36
Run the protein melt reactions	37
Load and run the reactions	37
Review the melt curves	38
Set up the analysis	40
Start the analysis software	40
Create a study	40
Experiment file guidelines	40
Set up the example ligand titration study	41
About the analysis settings	42
Review the analysis settings	42
Review the well results	42
About the melt curve plots	43
About the Boltzmann fit curve	44
Review the melt curves	44
About the analysis mode	45
Set the analysis mode	46
Review the regions of analysis (ROA)	46
Review the flags and T _m values	47
Review the Boltzmann fit	47
Example well results	48
Review the replicate results	48
About the T _m statistics	48
About the replicate results plot	48
Review the boltzmann T _m statistics	49
Review the positive hits	52
Example replicate results	54
Next steps	54

■ CHAPTER 5 Buffer screening studies	55
General guidelines	55
Experimental design	55
Replicates and controls	56
Create and set up an experiment file for the instrument run	56
Experiment file settings	56
Defining and assigning targets	57
Example targets	58
Example plate layout	58
Prepare the protein melt reactions	59
Protein melt reaction stability	59
Required materials	59
Prepare the protein melt reactions	59
Run the protein melt reactions	60
Load and run the reactions	60
Review the melt curves	61
Example melt curves	62
Set up the analysis	62
Setup guidelines	62
Create and set up the study	63
Review the analysis settings	66
Review the well results	66
Review the melt curves	67
Review the regions of analysis (ROA)	68
Review the flags and T _m values	69
Review the Boltzmann fit	70
Example well results	70
Review the replicate results	71
Review the T _m statistics	71
Review the positive hits	73

■ CHAPTER 6 Mutation screening studies	76
General guidelines	76
Experimental design	76
Replicates and controls	77
Create and set up an experiment file for the instrument run	77
Experiment file settings	77
Defining and assigning targets	78
Example targets	79
Example plate layout	79
Prepare the protein melt reactions	79
Protein melt reaction stability	79
Required materials	80
Prepare the protein melt reactions	80
Run the protein melt reactions	80
Load and run the reactions	81
Review the melt curves	81
Example melt curves	82
Set up the analysis	82
Setup guidelines	83
Create and set up the study	83
Review the analysis settings	85
Review the well results	85
Review the melt curves	86
Review the regions of analysis (ROA)	87
Review the flags and T _m values	88
Review the Boltzmann fit	89
Example well results	89
Review the replicate results	90
Review the T _m statistics	90
Review the positive hits	92

■	CHAPTER 7	Ligand screening studies	95
	General guidelines		95
	Experimental design		95
	Replicates and controls		96
	Create and set up an experiment file for the instrument run		96
	Experiment file settings		96
	Defining and assigning targets		97
	Example targets		98
	Example plate layout		98
	Prepare the protein melt reactions		98
	Protein melt reaction stability		98
	Required materials		98
	Prepare the protein melt reactions		99
	Run the protein melt reactions		99
	Load and run the reactions		100
	Review the melt curves		100
	Example melt curves		101
	Set up the analysis		101
	Setup guidelines		101
	Create and set up the study		102
	Review the analysis settings		104
	Review the well results		104
	Review the melt curves		105
	Review the regions of analysis (ROA)		106
	Review the flags and T _m values		107
	Review the Boltzmann fit		108
	Example well results		108
	Review the replicate results		109
	Review the T _m statistics		109
	Review the positive hits		111
■	CHAPTER 8	Troubleshooting	114
	Examples of observations		115
	High initial background signal		115
	Flat signal or decrease in signal		117
■	APPENDIX A	Supplemental information	118
	Hardware recommendations		118
	Boltzmann fitting method		119

■	APPENDIX B	Software warranty information	120
		Computer configuration	120
		Limited product warranty	120
		Limited warranty	120
		Warranty period effective date	120
		Warranty claims	121
		Warranty exceptions	121
		Warranty limitations	121
■	APPENDIX C	Ordering information	122
■	APPENDIX D	Glossary	123
■	APPENDIX E	Safety	126
		Chemical safety	127
		Biological hazard safety	128
■	APPENDIX F	Documentation and support	129
		Customer and technical support	129
		Limited product warranty	129



About this guide

IMPORTANT! Before using this product, read and understand the information the “Safety” appendix in this document.

Purpose

This guide is designed to help you quickly learn how to perform Protein Thermal Shift™ studies using Protein Thermal Shift™ reagents, Applied Biosystems™ Real-Time PCR Systems, and Protein Thermal Shift™ Software v1.4. This guide provides step-by-step procedures for the following tasks:

- How to perform an example ligand titration study using the Protein Thermal Shift™ Starter Kit and the example experiment files installed with the Protein Thermal Shift™ Software.
- How to perform and troubleshoot your buffer screening, mutation screening, and ligand screening studies using Protein Thermal Shift™ reagents and Protein Thermal Shift™ Software.

User attention words

Five user attention words may appear in this document. Each word implies a particular level of observation or action as described below:

Note: Provides information that may be of interest or help but is not critical to the use of the product.

IMPORTANT! Provides information that is necessary for proper instrument operation or accurate chemistry kit use.



CAUTION! Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.



WARNING! Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.



DANGER! Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury.



Getting started with a Protein Thermal Shift™ study

Perform a Protein Thermal Shift™ study to screen for buffers, mutations, or ligands that affect the thermal stability of the protein of interest. With Protein Thermal Shift™ studies, you can determine optimal protein storage conditions, perform functional characterization, or improve the success rates of protein purification and crystallization.

This chapter provides instructions for using the Protein Thermal Shift™ Starter Kit to perform an example ligand titration study. In the example ligand titration study, you titrate a ligand to determine the concentration that increases the thermal stability of the protein.

Workflow for a Protein Thermal Shift™ study

Create and set up an experiment file for the instrument run (page 15)

Prepare the reactions (page 36)

Run the protein melt reactions (page 37)

Set up the analysis (page 40)

Review the well results (page 42)

Review the replicate results (page 48)

Product information

Purpose of the starter kit

The Protein Thermal Shift™ Starter Kit includes buffer, dye, a control protein, and a control ligand.

After you follow the instructions in this chapter to learn how to perform a Protein Thermal Shift™ study, you can continue to use the starter kit components:

Use the Protein Thermal Shift™ Buffer and the Protein Thermal Shift™ Dye in your protein melt reactions. The buffer and dye included in the starter kit are identical to the buffer and dye in the Protein Thermal Shift™ Dye Kit.

Starter kit contents and storage

The Protein Thermal Shift™ Starter Kit (Cat. No. [4462263](#)) contains two boxes.

Box	Quantity	Component	Storage conditions
Protein Thermal Shift™ Dye Kit	Sufficient for 1000 reactions	Protein Thermal Shift™ Buffer	Room temperature (RT, 18°C to 25°C)
		Protein Thermal Shift™ Dye	
Protein Thermal Shift™ Starter Kit	Sufficient for 100 reactions	Protein Thermal Shift™ Control Ligand	–25°C to –15°C
		Protein Thermal Shift™ Control Protein	

Purpose of the dye kit

After you learn how to perform a Protein Thermal Shift™ study, you can use the Protein Thermal Shift™ Dye Kit. The Protein Thermal Shift™ Dye Kit contains buffer and dye to perform the Protein Thermal Shift™ reactions with your proteins of interest.

Dye kit contents and storage

The Protein Thermal Shift™ Dye Kit (Cat. No. [4461146](#)) contains one box.

Quantity	Component	Storage conditions
Sufficient for 1000 reactions	Protein Thermal Shift™ Buffer	Room temperature (RT, 18°C to 25°C)
	Protein Thermal Shift™ Dye	

Materials required but not included

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Catalog numbers that appear as links open the web pages for those products.

Item	Source
Applied Biosystems™ Real-Time PCR System	Contact your local sales office
Applied Biosystems™ Protein Thermal Shift™ Software v1.4	4466037
Applied Biosystems™ MicroAmp™ Optical 96-Well Reaction Plate with Barcode	4306737
Centrifuge with plate adapters	MLS
Microcentrifuge	MLS
Microcentrifuge tubes	MLS
Pipettors and pipette tips	MLS
Vortexer	MLS
Water	MLS

General guidelines

Protein considerations

Before you perform a Protein Thermal Shift™ study, consider whether the protein of interest is a suitable candidate for this method:

- Is the protein thermally stable, as determined by other methods? If the T_m for the protein is greater than 98°C, the protein melt reactions will produce melt curves with no distinct melt phase. You may observe flat fluorescence signals, a decrease in signal, or high signals, depending on the protein. You may need to use other methods to screen for conditions that increase thermal stability of the protein.

For thermally stable proteins, you may consider performing Protein Thermal Shift™ studies to screen for buffers, ligands, or mutations that *decrease* thermal stability of the protein.

- In the native state of a protein, are there external hydrophobic sites? If so, you may observe a high initial background signal. Conversely, if the protein does not contain sufficient hydrophobic residues, you may observe low fluorescence signals.
- Does the protein of interest comprise more than one domain or form oligomers? If so, you may observe a multi-state model of unfolding and multiple melt phases in the melt curves.
 - For proteins that comprise more than one domain, you may consider separate expression of just the domain with the active site.
 - For proteins that form oligomers, you can perform a buffer/additive screening study to identify conditions in which the protein unfolding follows a two-state model.

Experimental conditions

We recommend that you begin with the experimental conditions that we provide. If you do not observe clear, well-resolved melt phases in the melt curves and a reasonable rise in fluorescence relative to the NPC wells, you can optimize the experimental conditions according to your protein:

- **Protein:dye ratio:** If you observe high initial fluorescence or small transitional increase in your signal, then we recommend that you perform a titration study with the protein and Protein Thermal Shift™ Dye to identify the optimal protein:dye ratio.
- **Ramp speed and ramp rate:** Start with the Fast ramp speed and the recommended ramp rate, then decrease the ramp speed and/or ramp rate as needed to obtain well-resolved melt phases. In Protein Thermal Shift™ studies, the ramp speed affects protein unfolding and the resulting melt curves. With some proteins, we have observed low resolution of the melt phases in the melt curves with the Fast ramp speed and high ramp rates. For these proteins, we observed improved resolution of the melt phases when we used the Standard ramp speed and decreased ramp rates.
- **Thermal profile:** If you observe that the T_m is close to the beginning or close to the end of the melt curve, adjust the temperature range so that the T_m is in the middle of the temperature range and you are able to observe a complete protein melt.

Plate layout

Because of possible plate-to-plate variability, we recommend that you set up the plates so that each plate contains a reference group and an analysis group does not span multiple plates.

Replicates

For all Protein Thermal Shift™ studies, we recommend that you prepare at least 4 replicates of each protein melt reaction to ensure statistically significant results.

Controls

For Protein Thermal Shift™ studies, we recommend that you prepare controls:

- **No Protein Control (NPC):** Protein melt reactions that contain only buffer, water, and dye. NPC wells with high fluorescence signals and melt phases may indicate contamination in wells or protein melt reactions.
- **Ligand Only Control (LOC):** Protein melt reactions that contain only ligand, buffer, water, and dye. LOC wells with melt profiles distinct from NPC wells may indicate ligand-dye interactions. Ligands that bind the dye and affect the fluorescence levels may mask the presence or absence of protein-dye interactions.

Supported real-time PCR systems

Melt curve experiment files from the following Applied Biosystems™ Real-Time PCR Systems and system software are supported:

Applied Biosystems™ Real-Time PCR System	System software versions
QuantStudio™ 7 Pro Real-Time PCR System QuantStudio™ 6 Pro Real-Time PCR System	QuantStudio™ Design and Analysis Software v2 (v2.x)
QuantStudio™ 5 Real-Time PCR System QuantStudio™ 3 Real-Time PCR System QuantStudio™ 1 Real-Time PCR System	QuantStudio™ Design and Analysis Desktop Software (v1.x)
QuantStudio™ 7 Flex Real-Time PCR System QuantStudio™ 6 Flex Real-Time PCR System	QuantStudio™ Real-Time PCR Software (v1.x)
QuantStudio™ 12K Flex Real-Time PCR System	QuantStudio™ 12K Flex Software (v1.x)
ViiA™ 7 Real-Time PCR System	ViiA™ 7 Software v1.0 and v1.1
StepOne™ Real-Time PCR System and StepOnePlus™ Real-Time PCR System	StepOne™ Software v2.1 and v2.2
7500 Real-Time PCR System and 7500 Fast Real-Time PCR System	7500 Software v2.0.4 and v2.0.5

Note: Other system software versions for the real-time PCR systems listed may be accepted by the Protein Thermal Shift™ Software. By default, a warning message is displayed when you import an experiment file from a system software version that is not supported.



Create and set up an experiment file for the instrument run

This chapter describes how to set up an experiment file in the following software.





- QuantStudio™ Design and Analysis Software v2 (v2.x)
- QuantStudio™ Design and Analysis Desktop Software (v1.x)
- QuantStudio™ Real-Time PCR Software (v1.x)
- QuantStudio™ 12K Flex Software (v1.x)
- ViiA™ 7 Software
- StepOne™ Software
- 7500 Software

Set up a plate file for high resolution melt analysis in QuantStudio™ Design and Analysis Software v2 (v2.x)

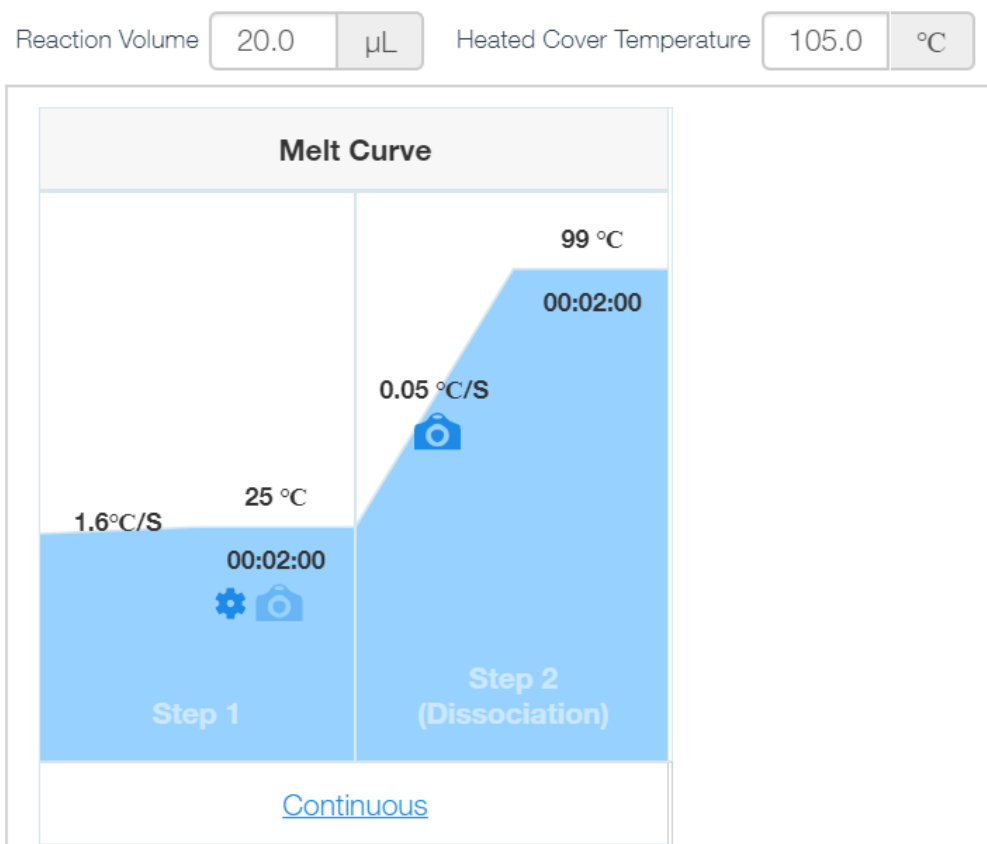
Define and assign targets so you can review the melt curves for the replicate groups in QuantStudio™ Design and Analysis Software v2 (v2.x).

This procedure is compatible with data files for the following real-time PCR instruments.

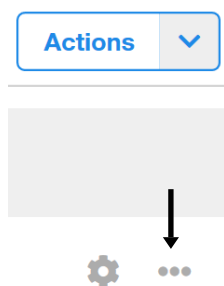
- QuantStudio™ 7 Pro Real-Time PCR System
- QuantStudio™ 6 Pro Real-Time PCR System

1. In the home screen, click  **Set Up Plate**.
2. Navigate to, then select a system template or plate file with a melt curve.
3. In a plate file, in the **Run Method** tab, adjust the run method.
 - a. Place the pointer over the **Hold** stage, then click .
 - b. Place the pointer over the **PCR** stage, then click .
 - c. In the **Melt Curve** stage, place the pointer over **Step 2**, then click .
 - d. Define the thermal profile.

Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	1.6°C/s	25.0	02:00
2	0.05°C/s	99.0	02:00



- Click ... (**Actions**) ▶ **Filter Settings**. In the **Filter settings** pane, click the **Melt Curve** tab, select the filters.



We recommend the following settings for optimal results.

Filter type	Selection	Example																																																	
Coupled	x2(520±10)—m2(558±11)	<div><div>Filter Settings</div><div><div>PCR</div><div>Melt Curve</div></div><table><thead><tr><th></th><th>m1(520±15)nm</th><th>m2(558±11)nm</th><th>m3(586±10)nm</th><th>m4(623±14)nm</th><th>m5(682±14)nm</th></tr></thead><tbody><tr><td>x1(470±15)nm</td><td><input type="checkbox"/></td><td></td><td></td><td></td><td></td></tr><tr><td>x2(520±10)nm</td><td></td><td><input checked="" type="checkbox"/></td><td></td><td></td><td></td></tr><tr><td>x3(550±11)nm</td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td></tr><tr><td>x4(580±10)nm</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td></tr><tr><td>x5(640±10)nm</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr></tbody></table><div><div>Cancel</div><div>Save</div></div></div>		m1(520±15)nm	m2(558±11)nm	m3(586±10)nm	m4(623±14)nm	m5(682±14)nm	x1(470±15)nm	<input type="checkbox"/>					x2(520±10)nm		<input checked="" type="checkbox"/>				x3(550±11)nm			<input type="checkbox"/>			x4(580±10)nm				<input type="checkbox"/>		x5(640±10)nm					<input type="checkbox"/>													
	m1(520±15)nm	m2(558±11)nm	m3(586±10)nm	m4(623±14)nm	m5(682±14)nm																																														
x1(470±15)nm	<input type="checkbox"/>																																																		
x2(520±10)nm		<input checked="" type="checkbox"/>																																																	
x3(550±11)nm			<input type="checkbox"/>																																																
x4(580±10)nm				<input type="checkbox"/>																																															
x5(640±10)nm					<input type="checkbox"/>																																														
Decoupled	x1(470±15)—m3(586±10)	<div><div>Filter Settings</div><div><div>PCR</div><div>Melt Curve</div></div><table><thead><tr><th></th><th>m1(520±15)nm</th><th>m2(558±11)nm</th><th>m3(586±10)nm</th><th>m4(623±14)nm</th><th>m5(682±14)nm</th><th>m6(711±12)nm</th></tr></thead><tbody><tr><td>x1(470±15)nm</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input checked="" type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x2(520±10)nm</td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x3(550±11)nm</td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x4(580±10)nm</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x5(640±10)nm</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x6(662±10)nm</td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr></tbody></table><div><div>Cancel</div><div>Save</div></div></div>		m1(520±15)nm	m2(558±11)nm	m3(586±10)nm	m4(623±14)nm	m5(682±14)nm	m6(711±12)nm	x1(470±15)nm	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x2(520±10)nm		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x3(550±11)nm			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x4(580±10)nm				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x5(640±10)nm					<input type="checkbox"/>	<input type="checkbox"/>	x6(662±10)nm						<input type="checkbox"/>
	m1(520±15)nm	m2(558±11)nm	m3(586±10)nm	m4(623±14)nm	m5(682±14)nm	m6(711±12)nm																																													
x1(470±15)nm	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																													
x2(520±10)nm		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																													
x3(550±11)nm			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																													
x4(580±10)nm				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																													
x5(640±10)nm					<input type="checkbox"/>	<input type="checkbox"/>																																													
x6(662±10)nm						<input type="checkbox"/>																																													

5. Add samples and targets, enter reagent information, and make the following selections.
 - a. In the upper-left corner of the **Plate Setup** tab, select **NONE** as the passive reference from the dropdown list.
 - b. In the **Targets** table, in the **Reporter** column, select ROX.

IMPORTANT! The software automatically assigns a task to the target in a well based on the sample type in that well. The **Positive Control** task is not compatible with Protein Thermal Shift™ Software; do not select **Positive Control** as the sample type.

You can use a single target and a single sample. Plate setup is completed later, in Protein Thermal Shift™ Software.





6. Review the plate file and send to the instrument run queue.

Set up a melt curve run in QuantStudio™ Design and Analysis Desktop Software (v1.x)

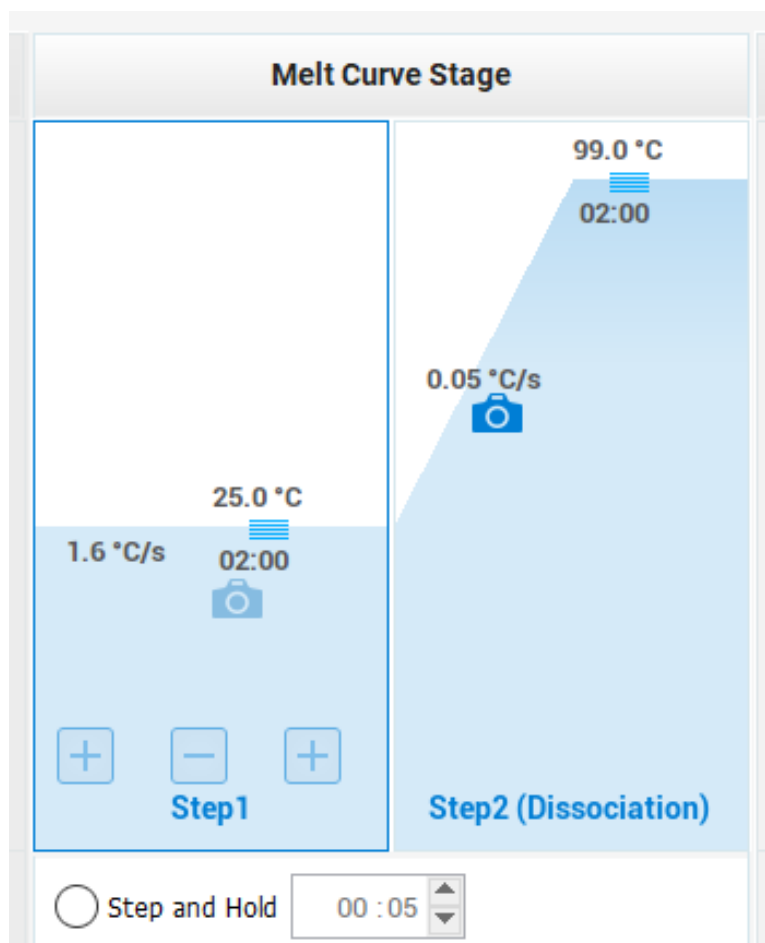
Define and assign targets so you can review the melt curves for the replicate groups in QuantStudio™ Design and Analysis Desktop Software (v1.x).

This procedure is compatible with data files for the following real-time PCR instruments.

- QuantStudio™ 5 Real-Time PCR System
- QuantStudio™ 3 Real-Time PCR System
- QuantStudio™ 1 Real-Time PCR System

1. In the  **Home** screen, create or open a template.
 - In the  **New Experiment** pane, click **Create New Experiment** to create a new template.
 - In the  **Open Existing Experiment** pane, click **Open** to select and open an existing template.
2. In the **Properties** tab, enter the template information.
 - a. In **Experiment type**, select **Melt Curve**.
 - b. In **Chemistry**, select **Other**.
 - c. Click **Next**.
3. In the **Method** tab, edit the default run method according to the run requirements.
 - a. Place the pointer over the **Step 2** stage, then click .
 - b. Define the thermal profile.

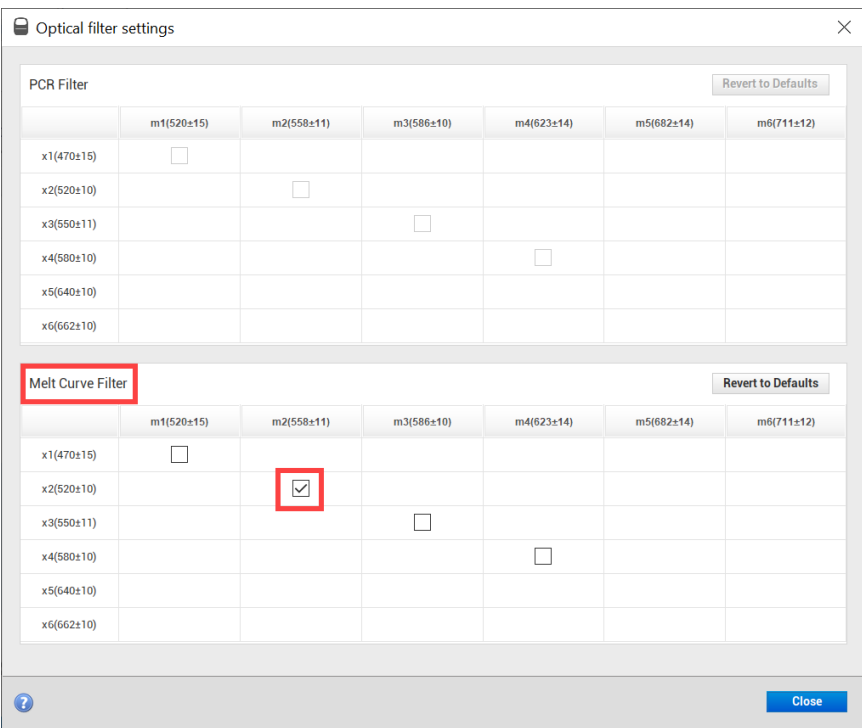
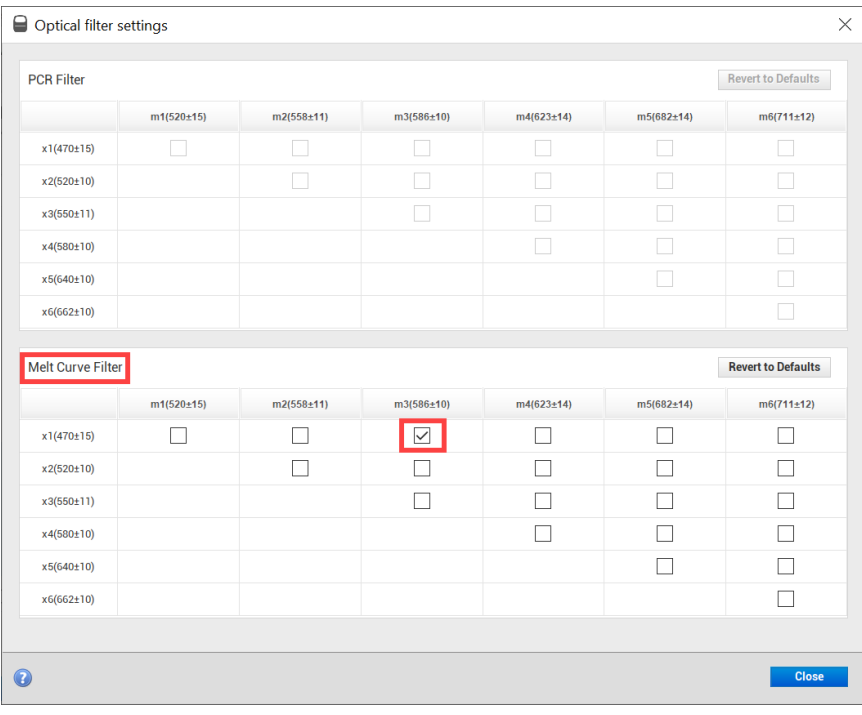
Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	1.6°C/s	25.0	02:00
2	0.05°C/s	99.0	02:00



- c. Select  **Action ▶ Optical filter settings.**



The excitation (x) and emission (m) wavelengths that correspond to each filter are shown on the screen.


- d. In the **Optical filter settings** pane, in the **Melt Curve Filter** section, select the filters.
We recommend the following settings for optimal results.

Filter type	Selection	Example
Coupled	x2(520±10) – m2(558±11)	 <p>The screenshot shows the 'Optical filter settings' dialog box. It has two main sections: 'PCR Filter' and 'Melt Curve Filter'. The 'Melt Curve Filter' section is highlighted with a red box. It contains a table with columns for filter types (m1 to m6) and rows for filter names (x1 to x6). The cell for x2(520±10) under the m2(558±11) column is checked, and this row is also highlighted with a red box.</p>
Decoupled	x1(470±15) – m3(586±10)	 <p>The screenshot shows the 'Optical filter settings' dialog box. It has two main sections: 'PCR Filter' and 'Melt Curve Filter'. The 'Melt Curve Filter' section is highlighted with a red box. It contains a table with columns for filter types (m1 to m6) and rows for filter names (x1 to x6). The cell for m3(586±10) under the x1(470±15) column is checked, and this cell is also highlighted with a red box.</p>

4. In the **Plate** tab (**Quick Setup**), assign plate attributes.
a. In the **Plate Attributes** pane, in the **Passive Reference** dropdown list, select **None**.

5. In the **Plate** tab (**Quick Setup**), define and assign well attributes.

- a. Select wells in the  **Plate Layout** or the  **Well Table**.
- b. Assign samples and targets to selected wells.
 - Enter new sample and target names in the text fields.
 - Select previously defined samples and targets from the dropdown lists.

Note: New sample or target names entered in the **Quick Setup** subtab are automatically populated with default values for **Reporter** (FAM) and **Quencher** (NFQ-MGB) and assigned a task ( **Unknown**). Edit these values in the **Advanced Setup** subtab.



6. (Optional) In the **Plate** tab (**Advanced Setup**), assign tasks.

- a. Select wells in the  **Plate Layout** or the  **Well Table**.

- b. In the **Targets** table, assign the targets.

You can use a single target. Plate setup is completed later, in Protein Thermal Shift™ Software.

1. Select the checkbox of a target, then select a task from the dropdown list.

Reaction type	Task
Unknown (test sample)	
No-template control	

2. In the Reporter dropdown list, select **ROX** for all wells that are filled.

Design an experiment in QuantStudio™ Real-Time PCR Software (v1.x)

Define and assign targets so that you can review the melt curves for the replicate groups in QuantStudio™ Real-Time PCR Software (v1.x). For real-time data collection, change the default analysis settings.

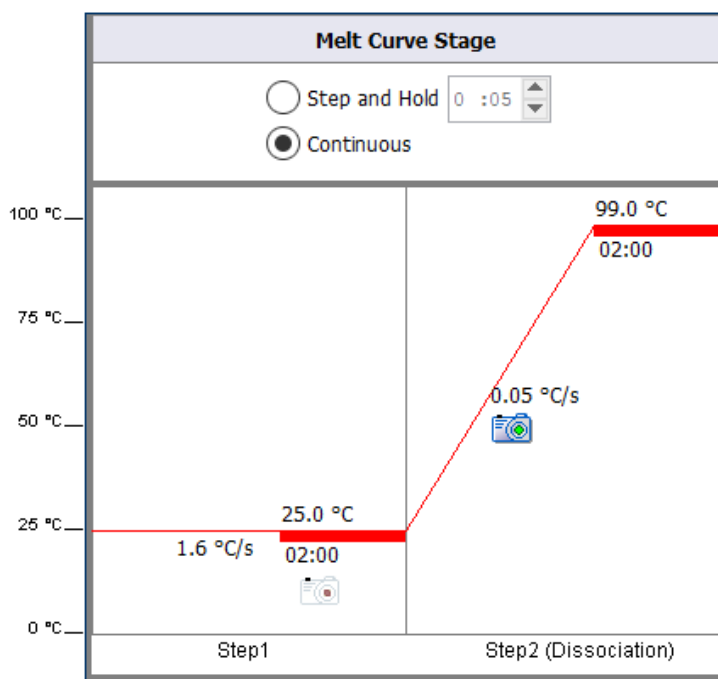
This procedure is compatible with data files for the following real-time PCR instruments.

- QuantStudio™ 7 Flex Real-Time PCR System
- QuantStudio™ 6 Flex Real-Time PCR System

1. Click **Tools ▶ Preferences**.
2. In the **Defaults tab**, select the **Show optical filters for run method** checkbox, then click **OK**.
3. Click **Experiment Setup**, then define the experiment properties. For the type of experiment, click **Melt Curve**.
4. Click **Setup ▶ Run Method** to set the thermal profile.
 - a. Select **Step 2**, then click **Delete Selected**.
 - b. Define the thermal profile.

We recommend the following conditions as a starting point. You can optimize the conditions.

Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	1.6°C/s	25.0	02:00
2	0.05°C/s	99.0	02:00



5. Click the **Optical Filters** tab. In the **Melt Curve Ramp Filter** section, select the filters.
We recommend the following settings for optimal results.

Filter type	Selection	Example																																																																																																		
Coupled	$x2(520\pm10) - m2(558\pm11)$	<div><div>Run Method</div><div>Reaction Volume per Well: <input type="text" value="20"/> μL</div><div><div>Graphical View</div><div>Optical Filters</div></div><div>PCR Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><thead><tr><th></th><th>m1(520\pm15)</th><th>m2(558\pm11)</th><th>m3(586\pm10)</th><th>m4(623\pm14)</th><th>m5(682\pm14)</th><th>m6(711\pm12)</th></tr></thead><tbody><tr><td>x1(470\pm15)</td><td><input type="checkbox"/></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x2(520\pm10)</td><td></td><td><input type="checkbox"/></td><td></td><td></td><td></td><td></td></tr><tr><td>x3(550\pm11)</td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td><td></td></tr><tr><td>x4(580\pm10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td></tr><tr><td>x5(640\pm10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td></tr><tr><td>x6(662\pm10)</td><td></td><td></td><td></td><td></td><td></td><td></td></tr></tbody></table><div>Melt Curve Ramp Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><thead><tr><th></th><th>m1(520\pm15)</th><th>m2(558\pm11)</th><th>m3(586\pm10)</th><th>m4(623\pm14)</th><th>m5(682\pm14)</th><th>m6(711\pm12)</th></tr></thead><tbody><tr><td>x1(470\pm15)</td><td><input type="checkbox"/></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x2(520\pm10)</td><td></td><td><input checked="" type="checkbox"/></td><td></td><td></td><td></td><td></td></tr><tr><td>x3(550\pm11)</td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td><td></td></tr><tr><td>x4(580\pm10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td></tr><tr><td>x5(640\pm10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td></tr><tr><td>x6(662\pm10)</td><td></td><td></td><td></td><td></td><td></td><td></td></tr></tbody></table></div>		m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)	x1(470 \pm 15)	<input type="checkbox"/>						x2(520 \pm 10)		<input type="checkbox"/>					x3(550 \pm 11)			<input type="checkbox"/>				x4(580 \pm 10)				<input type="checkbox"/>			x5(640 \pm 10)					<input type="checkbox"/>		x6(662 \pm 10)								m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)	x1(470 \pm 15)	<input type="checkbox"/>						x2(520 \pm 10)		<input checked="" type="checkbox"/>					x3(550 \pm 11)			<input type="checkbox"/>				x4(580 \pm 10)				<input type="checkbox"/>			x5(640 \pm 10)					<input type="checkbox"/>		x6(662 \pm 10)						
	m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)																																																																																														
x1(470 \pm 15)	<input type="checkbox"/>																																																																																																			
x2(520 \pm 10)		<input type="checkbox"/>																																																																																																		
x3(550 \pm 11)			<input type="checkbox"/>																																																																																																	
x4(580 \pm 10)				<input type="checkbox"/>																																																																																																
x5(640 \pm 10)					<input type="checkbox"/>																																																																																															
x6(662 \pm 10)																																																																																																				
	m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)																																																																																														
x1(470 \pm 15)	<input type="checkbox"/>																																																																																																			
x2(520 \pm 10)		<input checked="" type="checkbox"/>																																																																																																		
x3(550 \pm 11)			<input type="checkbox"/>																																																																																																	
x4(580 \pm 10)				<input type="checkbox"/>																																																																																																
x5(640 \pm 10)					<input type="checkbox"/>																																																																																															
x6(662 \pm 10)																																																																																																				
Decoupled	$x1(470\pm15) - m3(586\pm10)$	<div><div>Run Method</div><div>Reaction Volume per Well: <input type="text" value="20"/> μL</div><div><div>Graphical View</div><div>Optical Filters</div></div><div>PCR Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><thead><tr><th></th><th>m1(520\pm15)</th><th>m2(558\pm11)</th><th>m3(586\pm10)</th><th>m4(623\pm14)</th><th>m5(682\pm14)</th><th>m6(711\pm12)</th></tr></thead><tbody><tr><td>x1(470\pm15)</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x2(520\pm10)</td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x3(550\pm11)</td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x4(580\pm10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x5(640\pm10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x6(662\pm10)</td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr></tbody></table><div>Melt Curve Ramp Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><thead><tr><th></th><th>m1(520\pm15)</th><th>m2(558\pm11)</th><th>m3(586\pm10)</th><th>m4(623\pm14)</th><th>m5(682\pm14)</th><th>m6(711\pm12)</th></tr></thead><tbody><tr><td>x1(470\pm15)</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input checked="" type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x2(520\pm10)</td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x3(550\pm11)</td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x4(580\pm10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x5(640\pm10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x6(662\pm10)</td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr></tbody></table></div>		m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)	x1(470 \pm 15)	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x2(520 \pm 10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x3(550 \pm 11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x4(580 \pm 10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x5(640 \pm 10)					<input type="checkbox"/>	<input type="checkbox"/>	x6(662 \pm 10)						<input type="checkbox"/>		m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)	x1(470 \pm 15)	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x2(520 \pm 10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x3(550 \pm 11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x4(580 \pm 10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x5(640 \pm 10)					<input type="checkbox"/>	<input type="checkbox"/>	x6(662 \pm 10)						<input type="checkbox"/>
	m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)																																																																																														
x1(470 \pm 15)	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x2(520 \pm 10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x3(550 \pm 11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x4(580 \pm 10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x5(640 \pm 10)					<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x6(662 \pm 10)						<input type="checkbox"/>																																																																																														
	m1(520 \pm 15)	m2(558 \pm 11)	m3(586 \pm 10)	m4(623 \pm 14)	m5(682 \pm 14)	m6(711 \pm 12)																																																																																														
x1(470 \pm 15)	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x2(520 \pm 10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x3(550 \pm 11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x4(580 \pm 10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x5(640 \pm 10)					<input type="checkbox"/>	<input type="checkbox"/>																																																																																														
x6(662 \pm 10)						<input type="checkbox"/>																																																																																														

6. Click **Define**, then make the following selections.
 - a. In the **Passive Reference** list, select **none**.
 - b. In the **Targets** section, in the **Reporter** list, select **ROX** for each target.
 - c. Create one sample only.
The complete setup is done in Protein Thermal Shift™ Software.
7. Assign samples and ensure that all filled wells are assigned.
8. Save the experiment.

Design an experiment in QuantStudio™ 12K Flex Software

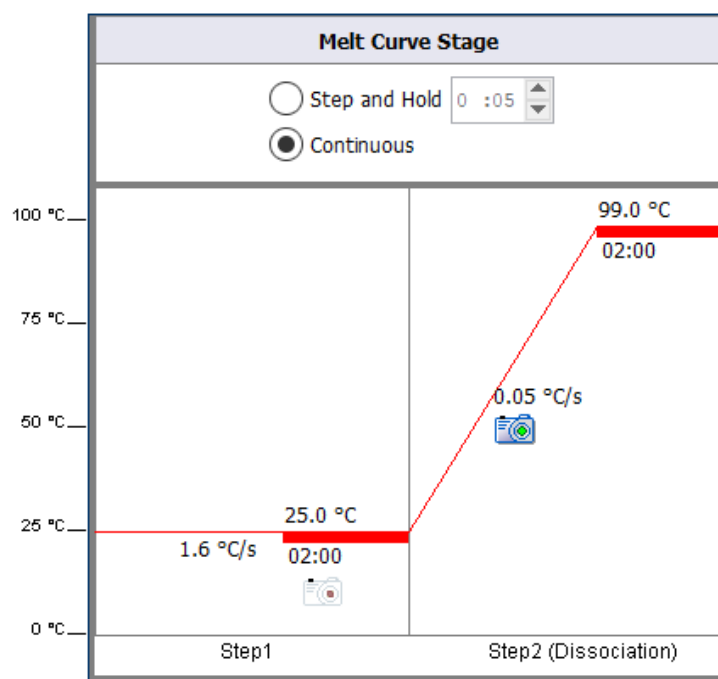
Define and assign targets so that you can review the melt curves for the replicate groups in QuantStudio™ 12K Flex Software. For real-time data collection, change the default analysis settings.

This procedure is compatible with data files for QuantStudio™ 12K Flex Real-Time PCR System.

1. Click **Tools ▶ Preferences**.
2. In the **Defaults tab**, select the **Show optical filters for run method** checkbox, then click **OK**.
3. Click **Experiment Setup**, then define the experiment properties. For the type of experiment, click **Melt Curve**.
4. Click **Setup ▶ Run Method** to set the thermal profile.
 - a. Select **Step 2**, then click **Delete Selected**.
 - b. Define the thermal profile.

We recommend the following conditions as a starting point. You can optimize the conditions.

Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	1.6°C/s	25.0	02:00
2	0.05°C/s	99.0	02:00



5. Click the **Optical Filters** tab. In the **Melt Curve Ramp Filter** section, select the filters.
We recommend the following settings for optimal results.

Filter type	Selection	Example																																																																																																																
Coupled	x2(520±10)—m2(558±11)	<div><div>Run Method</div><div>Reaction Volume per Well: 20 μL</div><div><div>Graphical View</div><div>Optical Filters</div></div><div>PCR Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><tr><td></td><td>m1(520±15)</td><td>m2(558±11)</td><td>m3(586±10)</td><td>m4(623±14)</td><td>m5(682±14)</td><td>m6(711±12)</td></tr><tr><td>Excitation Filter</td><td></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x1(470±15)</td><td><input type="checkbox"/></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x2(520±10)</td><td></td><td><input type="checkbox"/></td><td></td><td></td><td></td><td></td></tr><tr><td>x3(550±11)</td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td><td></td></tr><tr><td>x4(580±10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td></tr><tr><td>x5(640±10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td></tr><tr><td>x6(662±10)</td><td></td><td></td><td></td><td></td><td></td><td></td></tr></table><div>Melt Curve Ramp Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><tr><td></td><td>m1(520±15)</td><td>m2(558±11)</td><td>m3(586±10)</td><td>m4(623±14)</td><td>m5(682±14)</td><td>m6(711±12)</td></tr><tr><td>Excitation Filter</td><td></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x1(470±15)</td><td><input type="checkbox"/></td><td><input checked="" type="checkbox"/></td><td></td><td></td><td></td><td></td></tr><tr><td>x2(520±10)</td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td><td></td></tr><tr><td>x3(550±11)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td><td></td></tr><tr><td>x4(580±10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td></td></tr><tr><td>x5(640±10)</td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr><tr><td>x6(662±10)</td><td></td><td></td><td></td><td></td><td></td><td></td></tr></table></div>		m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)	Excitation Filter							x1(470±15)	<input type="checkbox"/>						x2(520±10)		<input type="checkbox"/>					x3(550±11)			<input type="checkbox"/>				x4(580±10)				<input type="checkbox"/>			x5(640±10)					<input type="checkbox"/>		x6(662±10)								m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)	Excitation Filter							x1(470±15)	<input type="checkbox"/>	<input checked="" type="checkbox"/>					x2(520±10)			<input type="checkbox"/>				x3(550±11)				<input type="checkbox"/>			x4(580±10)					<input type="checkbox"/>		x5(640±10)						<input type="checkbox"/>	x6(662±10)						
	m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)																																																																																																												
Excitation Filter																																																																																																																		
x1(470±15)	<input type="checkbox"/>																																																																																																																	
x2(520±10)		<input type="checkbox"/>																																																																																																																
x3(550±11)			<input type="checkbox"/>																																																																																																															
x4(580±10)				<input type="checkbox"/>																																																																																																														
x5(640±10)					<input type="checkbox"/>																																																																																																													
x6(662±10)																																																																																																																		
	m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)																																																																																																												
Excitation Filter																																																																																																																		
x1(470±15)	<input type="checkbox"/>	<input checked="" type="checkbox"/>																																																																																																																
x2(520±10)			<input type="checkbox"/>																																																																																																															
x3(550±11)				<input type="checkbox"/>																																																																																																														
x4(580±10)					<input type="checkbox"/>																																																																																																													
x5(640±10)						<input type="checkbox"/>																																																																																																												
x6(662±10)																																																																																																																		
Decoupled	x1(470±15)—m3(586±10)	<div><div>Run Method</div><div>Reaction Volume per Well: 20 μL</div><div><div>Graphical View</div><div>Optical Filters</div></div><div>PCR Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><tr><td></td><td>m1(520±15)</td><td>m2(558±11)</td><td>m3(586±10)</td><td>m4(623±14)</td><td>m5(682±14)</td><td>m6(711±12)</td></tr><tr><td>Excitation Filter</td><td></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x1(470±15)</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x2(520±10)</td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x3(550±11)</td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x4(580±10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x5(640±10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x6(662±10)</td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr></table><div>Melt Curve Ramp Filter</div><div><div>Load</div><div>Save</div><div>Revert to Defaults</div></div><div>Emission Filter</div><table><tr><td></td><td>m1(520±15)</td><td>m2(558±11)</td><td>m3(586±10)</td><td>m4(623±14)</td><td>m5(682±14)</td><td>m6(711±12)</td></tr><tr><td>Excitation Filter</td><td></td><td></td><td></td><td></td><td></td><td></td></tr><tr><td>x1(470±15)</td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input checked="" type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x2(520±10)</td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x3(550±11)</td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x4(580±10)</td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x5(640±10)</td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td><td><input type="checkbox"/></td></tr><tr><td>x6(662±10)</td><td></td><td></td><td></td><td></td><td></td><td><input type="checkbox"/></td></tr></table></div>		m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)	Excitation Filter							x1(470±15)	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x2(520±10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x3(550±11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x4(580±10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x5(640±10)					<input type="checkbox"/>	<input type="checkbox"/>	x6(662±10)						<input type="checkbox"/>		m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)	Excitation Filter							x1(470±15)	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x2(520±10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x3(550±11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x4(580±10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	x5(640±10)					<input type="checkbox"/>	<input type="checkbox"/>	x6(662±10)						<input type="checkbox"/>
	m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)																																																																																																												
Excitation Filter																																																																																																																		
x1(470±15)	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x2(520±10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x3(550±11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x4(580±10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x5(640±10)					<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x6(662±10)						<input type="checkbox"/>																																																																																																												
	m1(520±15)	m2(558±11)	m3(586±10)	m4(623±14)	m5(682±14)	m6(711±12)																																																																																																												
Excitation Filter																																																																																																																		
x1(470±15)	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x2(520±10)		<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x3(550±11)			<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x4(580±10)				<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x5(640±10)					<input type="checkbox"/>	<input type="checkbox"/>																																																																																																												
x6(662±10)						<input type="checkbox"/>																																																																																																												

6. Click **Define**, then make the following selections.
 - a. In the **Passive Reference** list, select **none**.
 - b. In the **Targets** section, in the **Reporter** list, select **ROX** for each target.
 - c. Create one sample only.
The complete setup is done in Protein Thermal Shift™ Software.
7. Assign samples and ensure that all filled wells are assigned.
8. Save the experiment.

Create and set up an experiment file for the ViiA™ 7 Real-Time PCR System

Define and assign targets so you can review the melt curves for the replicate groups in the ViiA™ 7 Software.

1. Select **Start ▶ All Programs ▶ Applied Biosystems™ ▶ ViiA™ 7 Software ▶ ViiA™ 7 Software v1.1.**

Note: If this is your first time starting the software, read and accept the license agreement.

2. Show the optical filters.

The **Optical Filters** tab is not visible by default. Enable the tab so that you can change the filter settings.

- a. Click **Tools ▶ Preferences.**
- b. In the **Defaults** tab, select the **Show optical filters for run method** checkbox, then click **OK.**

3. In the Home screen of the ViiA™ 7 Software, click **Experiment Setup.**






4. Complete the Experiment Properties screen:

Field	Entry
Experiment Name	Enter a unique Experiment Name using up to 100 letters and/or numbers. Note: If you plan to start the run using the instrument touchscreen, do not enter more than 32 characters for the Experiment Name and do not include spaces in the name.
Block type	Select the type of block that you are using: 384-Well Block or 96-Well Block (0.2mL)
Experiment type	Melt Curve
Reagent type	Other
Ramp speed	Fast

5. In the Define screen, define the targets (conditions) and dyes:

- a. In the Targets pane, enter the target name for each condition, select **ROX™** for the reporter, and select **None** for the Quencher.

For the example ligand titration study:

Target Name	Reporter	Quencher	Color
ProteinA +0 mM L	ROX	None	
ProteinA +0.1 mM L	ROX	None	
ProteinA +1 mM L	ROX	None	
LOC	ROX	None	
NPC	ROX	None	

- b. For the Passive Reference, select **None.**

6. In the Assign screen, assign the target and task to wells:
 - a. For each target, select the wells that contain the control protein, select the **Assign** checkbox for Control Protein, then select **U** for the Task.
 - b. (Optional) For no protein controls (NPC), select the NPC wells, select the **Assign** checkbox for Control Protein, then select **N** for the Task.
7. Ensure that the well assignments correspond exactly with the arrangement of reactions in the reaction plate.

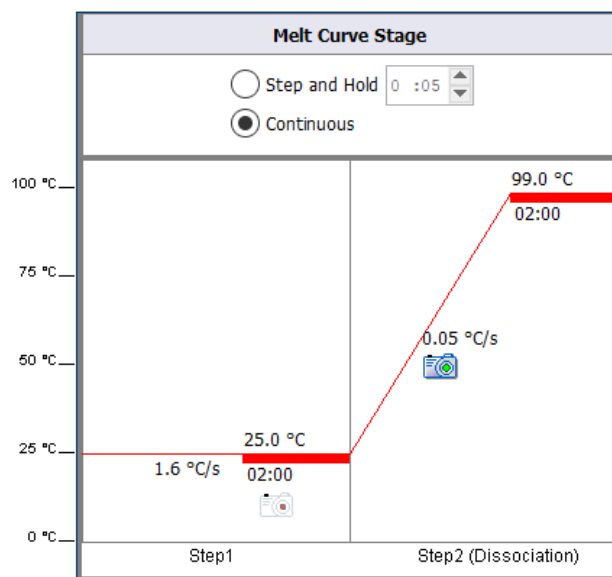
For the example ligation titration study:

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U
B	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U
C	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U	U
D	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc	U loc
E	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC

IMPORTANT! Setup errors can result in an incorrect grouping of replicates.

8. Complete the Run Method screen to define the melt curve:
 - a. For the Reaction Volume Per Well, enter **20** (μL).
 - b. For the ramp mode, select **Continuous**.
 - c. Define the thermal profile:

Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	1.6°C/s	25.0	02:00
2	0.05°C/s	99.0	02:00



9. Select the optical filters for the melt curve.

- a. In the **Run Method** screen, select the **Optical Filters** tab.

Note: If you do not see the Optical Filters tab, select **Tools ▶ Preferences**, then select the **Show optical filters for run method** checkbox in the Defaults tab.

- b. In the **Melt Curve Filter** section, make one of the following selections for the Excitation Filter-Emission Filter.

We recommend the following settings for optimal results.

- For coupled filters, select the checkbox in the row of **x2(520±10)** and the column of **m2(558±11)**.
- For decoupled filters, select the checkbox in the row of **x1(470±15)** and the column of **m3(586±10)**.


10. Select **File ▶ Save**, then enter a file name and select a location for the experiment file.

Create and set up an experiment file for the StepOne™ Real-Time PCR System or StepOnePlus™ Real-Time PCR System

Define and assign targets so you can review the melt curves for the replicate groups in the StepOne™ Software.

1. In the Home screen of StepOne™ Software, click **Advanced Setup**.

Note:

- If you are setting up the experiment file on a computer that is not connected to the instrument, click **Continue without Connection** in the Instrument Connection Failed dialog box.
- If you do not see the Advanced Setup button, click  below the Design Wizard button.






2. Complete the Experiment Properties screen:

Field	Entry
Experiment Name	Enter a unique Experiment Name using up to 50 letters and/or numbers.
Instrument type	Select the type of instrument that you are using: StepOnePlus™ Instrument (96 Wells) or StepOne™ Instrument (48 Wells)
Experiment type	Melt Curve
Reagent type	Other
Ramp speed	Fast

3. In the Plate Setup > Define Targets and Samples screen, define the targets (conditions):

In the Targets pane, enter the target name for each condition, select **ROX™** for the reporter, and select **None** for the Quencher.

For the example ligand titration study:

Target Name	Reporter	Quencher	Color
Protein + 0 mM L	ROX	None	
Protein+0.1mM L	ROX	None	
Protein +1mM L	ROX	None	
LOC	ROX	None	
NPC	ROX	None	

4. Define the well contents in the Plate Setup > Assign Targets and Samples screen:

- a. In the Plate Setup screen, select the **Assign Targets and Samples** tab.
- b. For each target (conditions), select the wells with those conditions, select the **Assign** checkbox for the target, then select **U** for the Task.

- c. (Optional) For no protein controls (NPC), select the NPC wells, select the **Assign** checkbox for the **NPC** target, then select **N** for the Task.
 - d. For the passive reference, select **None**.
- For the example ligand titration study (column 12 is empty):

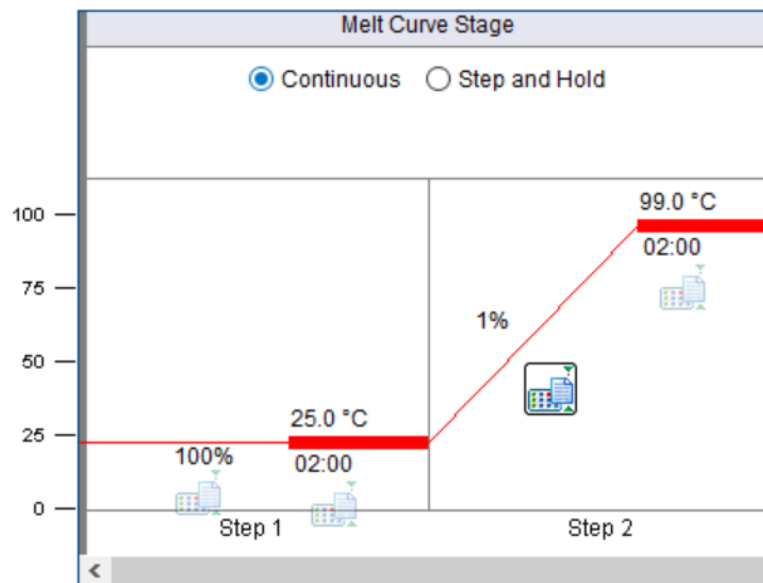
	1	2	3	4	5	6	7	8	9	10	11
A	U	U	U	U	U	U	U	U	U	U	U
B	U	U	U	U	U	U	U	U	U	U	U
C	U	U	U	U	U	U	U	U	U	U	U
D	U	U	U	U	U	U	U	U	U	U	U
E	U	U	U	U	U	U	U	U	U	U	U
F	U	U	U	U	U	U	U	U	U	U	U
G	U	U	U	U	U	U	U	U	U	U	U
H	U	U	U	U	U	U	U	U	U	U	U

5. Make sure that the well assignments correspond exactly with the arrangement of reactions in the reaction plate.

IMPORTANT! Setup errors may result in an incorrect grouping of replicates.

6. Complete the Run Method screen to define the melt curve:
 - a. For the Reaction Volume Per Well, enter **20** (μL).
 - b. For the ramp mode, select **Continuous**.
 - c. Define the thermal profile:

Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	100%	25.0	02:00
2	1%	99.0	02:00




7. Select **File ▶ Save**, then enter a file name and select a location for the experiment file.

Create and set up an experiment file for the 7500 Real-Time PCR System or 7500 Fast Real-Time PCR System

Define and assign targets so you can review the melt curves for the replicate groups in the 7500 Software.

1. Select **Start** ► **All Programs** ► **Applied Biosystems™** ► **7500 Software v2.0.5** to start the instrument software, then log into the software.
2. In the Home screen of the 7500 Software, click **Advanced Setup**.






Note: If you do not see the Advanced Setup button, click  below the Design Wizard button.

3. Complete the Experiment Properties screen:

Field	Entry
Experiment Name	Enter a unique Experiment Name using up to 100 letters and/or numbers.
Instrument type	7500 Fast (96 Wells) or 7500 (96 Wells)
Experiment type	Melt Curve
Reagent type	Other
Ramp speed	Standard

4. In the Plate Setup > Define Targets and Samples screen, define the targets (conditions):
In the Targets pane, enter the target name for each condition, select **ROX™** for the reporter, and select **None** for the Quencher.

For the example ligand titration study:

Target Name	Reporter	Quencher	Color
T4 lig	ROX	None	
T4lig+ 0.1mM L	ROX	None	
T4lig+ 1mM L	ROX	None	
NPC	ROX	None	
LOC	ROX	None	

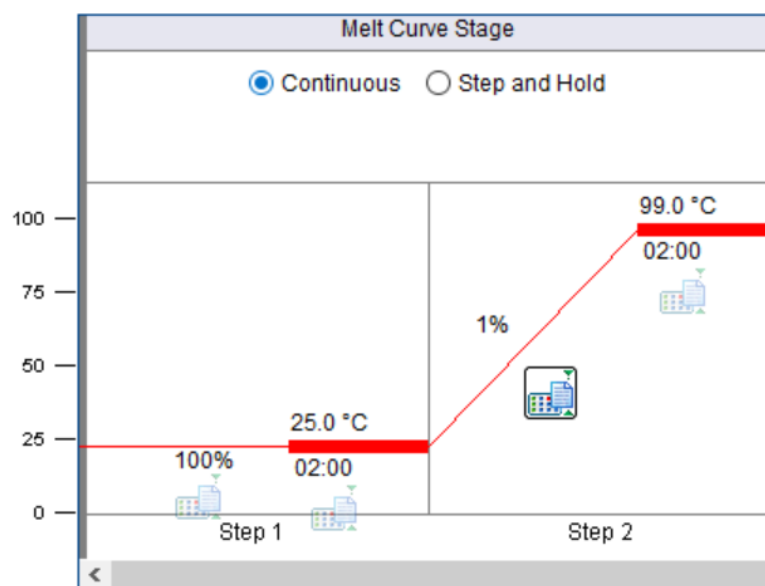
5. Define the well contents in the Plate Setup > Assign Targets and Samples screen:
 - a. In the Plate Setup screen, click the **Assign Targets and Samples** tab.
 - b. For each target (conditions), select the wells with those conditions, select the **Assign** checkbox for the target, then select **U** for the Task.
 - c. (Optional) For no protein controls (NPC), select the NPC wells, select the **Assign** checkbox for the **NPC** target, then select **N** for the Task.
 - d. For the passive reference, select **None**.

6. Make sure that the well assignments correspond exactly with the arrangement of reactions in the reaction plate.

IMPORTANT! Setup errors may result in an incorrect grouping of replicates.

7. Complete the Run Method screen to define the melt curve:
 - a. For the Reaction Volume Per Well, enter **20** (μL).
 - b. For the ramp mode, select **Continuous**.
 - c. Define the thermal profile:

Step	Ramp rate	Temp (°C)	Time (mm:ss)
1	100%	25.0	02:00
2	1%	99.0	02:00



8. Select **File ▶ Save**, then enter a file name and select a location for the experiment file.



Prepare and run the reactions, then review results

About the example study protein melt reactions

For the example ligand titration study, prepare 5 sets of protein melt reactions using the control protein and control ligand in the Protein Thermal Shift™ Starter Kit:

- Control protein and 0.0 mM control ligand
- Control protein and 0.1 mM control ligand
- Control protein and 1.0 mM control ligand
- No protein and 0.0 mM control ligand (NPC)
- No protein and 1.0 mM control ligand (LOC)

Required materials

Required materials for protein melt reactions:

- Protein Thermal Shift™ Dye (1000X)
- Protein Thermal Shift™ Buffer
- Water
- Protein Thermal Shift™ Control Protein
- Protein Thermal Shift™ Control Ligand
- MicroAmp™ Optical Reaction Plate appropriate for your real-time PCR instrument
- MicroAmp™ Optical Adhesive Film appropriate for your reaction plate

Prepare the protein melt reactions

We recommend that you prepare at least four replicates of each reaction.

1. Prepare a fresh dilution of Protein Thermal Shift™ Dye (1000X) to 8X.
2. Place the appropriate reaction plate or tubes on ice, then prepare the protein melt reactions:
 - Make sure that the arrangement of reactions in the reaction plate corresponds exactly with the well assignments in the experiment file.

IMPORTANT! Setup errors may result in an incorrect grouping of the data and incorrect T_m statistics.

- Add reaction components to the plate in the order listed.

Component	Volume
Protein Thermal Shift™ Buffer	5.0 µL
Water + 2.0 µL Protein Thermal Shift™ Control Protein + Protein Thermal Shift™ Control Ligand (0 mM, 0.1 mM, or 1 mM final concentration)	12.5 µL
Diluted Protein Thermal Shift™ Dye (8X)	2.5 µL
Total volume for each control reaction	20.0 µL

3. Pipet each reaction up and down 10 times to mix well.
4. Seal the plate with MicroAmp™ Optical Adhesive Film, spin it at 1000 rpm for 1 minute, then place it on ice.

Run the protein melt reactions

Load and run the protein melt reactions on a supported Applied Biosystems™ Real-Time PCR System, then analyze and save the experiment file before importing the experiment file into the Protein Thermal Shift™ Software.

IMPORTANT! Keep the protein melt reactions on ice until you load the instrument.

Load and run the reactions

For instructions about how to transfer experiment files to the instrument computer or instrument, how to operate the instrument, how to start the run from an instrument touchscreen, and how to monitor the run, see the instrument user guide.

IMPORTANT! Do not set the temperature below room temperature when you load and start the instrument. Damage to the instrument can occur if the temperature is below room temperature during a run.

1. Transfer the experiment file that you created for the run to the computer that is connected to the instrument, if needed.
2. In the Home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file you created for the run.
3. Load the reaction plate into the instrument.
4. In the Real-Time PCR System Software, click **Run** in the navigation pane, then click **START RUN**.

Review the melt curves

Using the Real-Time PCR System Software, open the experiment file from the completed instrument run, analyze and save the experiment file, then review the melt curves.

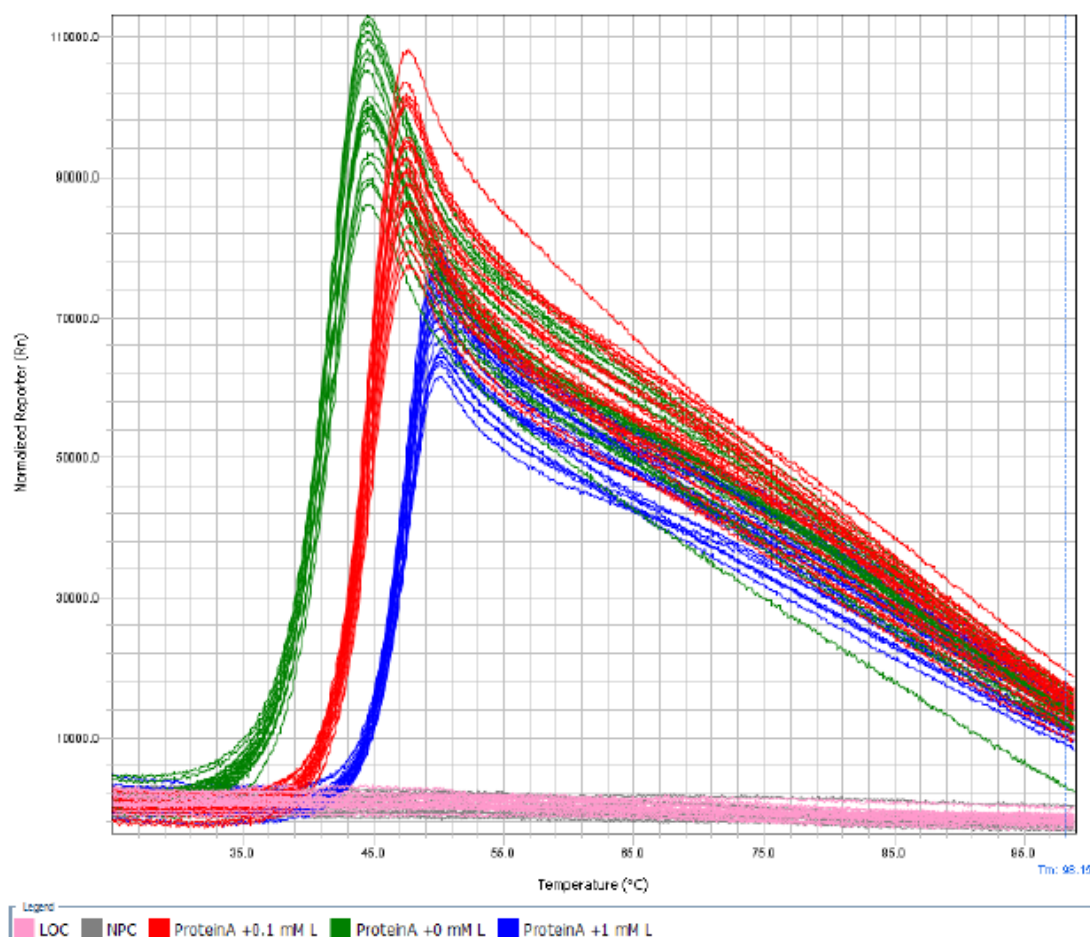
Note: You must analyze and save the experiment file in the Real-Time PCR System Software before you can import it into the Protein Thermal Shift™ Software.

Some common troubleshooting causes are provided here. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

1. In the Home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file from the instrument run.
2. View the melt curves:

Real-Time PCR System Software	View the melt curve
QuantStudio™ Design and Analysis Software v2 (v2.x)	<ol style="list-style-type: none"> 1. In the Quality Check tab, in the plot pane, select Melt Curve Plot from the dropdown list. 2. In the plot pane, click ⚙️ (Settings), then make the following selections. <ul style="list-style-type: none"> • Color By: Target • Plot Type: Normalized.
QuantStudio™ Design and Analysis Desktop Software (v1.x)	<ol style="list-style-type: none"> 1. In the Results tab, select Melt Curve Plot from the dropdown list. 2. Click 👁️ to configure the plot with the following selections. <ul style="list-style-type: none"> • Plot Type: Normalized • Color: Target. • Target: All or a target of interest
QuantStudio™ Real-Time PCR Software (v1.x) and QuantStudio™ 12K Flex Software (v1.x)	<ol style="list-style-type: none"> 1. Click Analysis ▶ Melt Curve Plot. 2. In the plot pane, make the following selections. <ul style="list-style-type: none"> • Color: Target • Plot: Normalized Reporter.
ViiA™ 7 Software	Click Analysis ▶ Melt Curve Plot in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.
StepOne™ Software or 7500 Software	Click Analysis ▶ Melt Curve in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.

Melt curves for the ligand screening example file from the ViiA™ 7 Real-Time PCR System:



3. Review the melt curves:

- Do you see fluorescence signals in all of the sample wells?
No fluorescence signals in the sample wells may indicate missing dye or protein or an instrument problem.
- Do you see flat fluorescence levels in the NPC wells?
High fluorescence levels in the NPC wells may indicate protein contamination in the wells or protein melt reactions; or it may indicate that the dye interacts with a component in the buffer.
- Do you see flat fluorescence levels in the LOC wells?
With the Protein Thermal Shift™ Control Ligand, high fluorescence levels in the LOC wells but not in the NPC wells may indicate protein contamination in the control ligand.
With your own samples, high fluorescence levels in the LOC wells but not in the NPC wells may indicate protein contamination in the control ligand or ligand-dye interactions.
- Do the replicates have similar melt curves?

4. Save, then close the experiment file.

Note: The melt curves in the real-time PCR software may not exactly match the melt curves in the Protein Thermal Shift™ Software. When the experiment files are imported into the Protein Thermal Shift™ Software, the Protein Thermal Shift™ Software reduces the noise in the fluorescence data.


Set up the analysis

This section provides instructions for setting up the Protein Thermal Shift™ study using Protein Thermal Shift™ Software v1.4 and the example ligand screening files that are installed with the software.

Start the analysis software

Select **Start ▶ All Programs ▶ Applied Biosystems™ ▶ Protein Thermal Shift™ Software v1.4**.

Create a study

1. In the Home screen, click  **Create Study** in the toolbar.
2. In the Setup > Properties screen, enter a unique name for the study.

Note: The Study Name cannot be more than 100 characters and cannot contain these characters: / \ * " ? < > | . ,

3. For the Instrument, select the instrument that you used to run the protein melt reactions.

IMPORTANT! The instrument selection must match the instrument type that you used to run the protein melt reactions and generate the experiment files.

Experiment file guidelines

Guidelines for the experiment files that you import into the study:


- The experiment file must contain analyzed and saved melt curve data from a complete melt curve run.
- The instrument type in the experiment file must match the instrument type selected for the study.

Note: The data needed for the study are extracted from the EDS file. The Protein Thermal Shift™ Software does not create, modify, save, or export EDS files.


Set up the example ligand titration study

First import the example experiment file (EDS), then set up the plate using the example plate template file (CSV).

The example files are located in <...>Program Files\Applied Biosystems\Protein Thermal Shift Software\examples, where <...> is the installation drive.

1. Go to the Setup > Experiment Files screen, click  **Import**, then select the ligand screening example file (EDS) for the instrument type that you selected for the study:
 - Ligand_Screening_Example_ViiA7.eds
 - Ligand_Screening_Example_StepOnePlus.eds
 - Ligand_Screening_Example_7500.eds

Note: The experiment file name and data are imported. Plate setup information is not imported.

2. Click  **Load Plate Template**, then select the ligand screening plate template file (CSV) for the instrument type that you selected for the study:
 - Ligand_Screening_Example_ViiA7.csv
 - Ligand_Screening_Example_StepOnePlus.csv
 - Ligand_Screening_Example_7500.csv



The conditions, task, and analysis group are assigned to the wells that contained protein melt reactions.

3. Make sure that the condition assignments correspond exactly with the contents of the reaction plate.

IMPORTANT! Setup errors may result in an incorrect grouping of the data and incorrect T_m statistics.

4. Select the replicate group to use as the reference:
 - a. Select the wells that contain 0.1 mM ligand:

Instrument type	Select wells
ViiA™ 7 Real-Time PCR System	Click the B row header to select row B (wells B1-B24).
StepOnePlus™ Real-Time PCR System or 7500 Fast Real-Time PCR System	Click-drag the 4 column header to select columns 46 (wells A4-A6, B4-B6, C4-C6, D4-D6, E4-E6, F4-F6, G4-G6, H4-H6).

- b. Click  **Assign**, then select **Reference** as the Task.
5. Click  **Save** in the toolbar to save and analyze the study.


About the analysis settings

With the analysis settings, you can adjust how positive hits are determined and specify settings for applying flags. If you generate a study template from the study file, the analysis settings for the study are saved in the study template.

Although the flags and the positive hits can be useful ways to quickly scan the results, we recommend that you review the values in the tables, review the plots, and review the replicates carefully.

Review the analysis settings

Review and edit the analysis settings to optimize the analysis for your study.

1. Click  **Analysis Settings** in the toolbar.
2. On the Positive Hit tab, specify the ΔT_m -Boltzmann and ΔT_m -Derivative values to indicate a positive hit:
 - Select > to identify conditions that increase protein thermal stability, or select < to identify conditions that decrease protein thermal stability.
 - Enter the number of degrees (°C) of T_m shift relative to the reference to indicate a positive hit.
3. On the Flags tab, specify settings for applying flags:
 - a. Select the flags to use in the analysis.
 - b. For the High Background, High NPC, Low Signal, and Poor Fit flags, specify the condition and threshold for applying the flag.
4. Apply the analysis settings and analyze:
 - Click **Apply** to apply the analysis settings and reanalyze while keeping the Analysis Settings dialog box open.
or
 - Click **OK** to apply the analysis settings, reanalyze, and close the Analysis Settings.

Review the well results

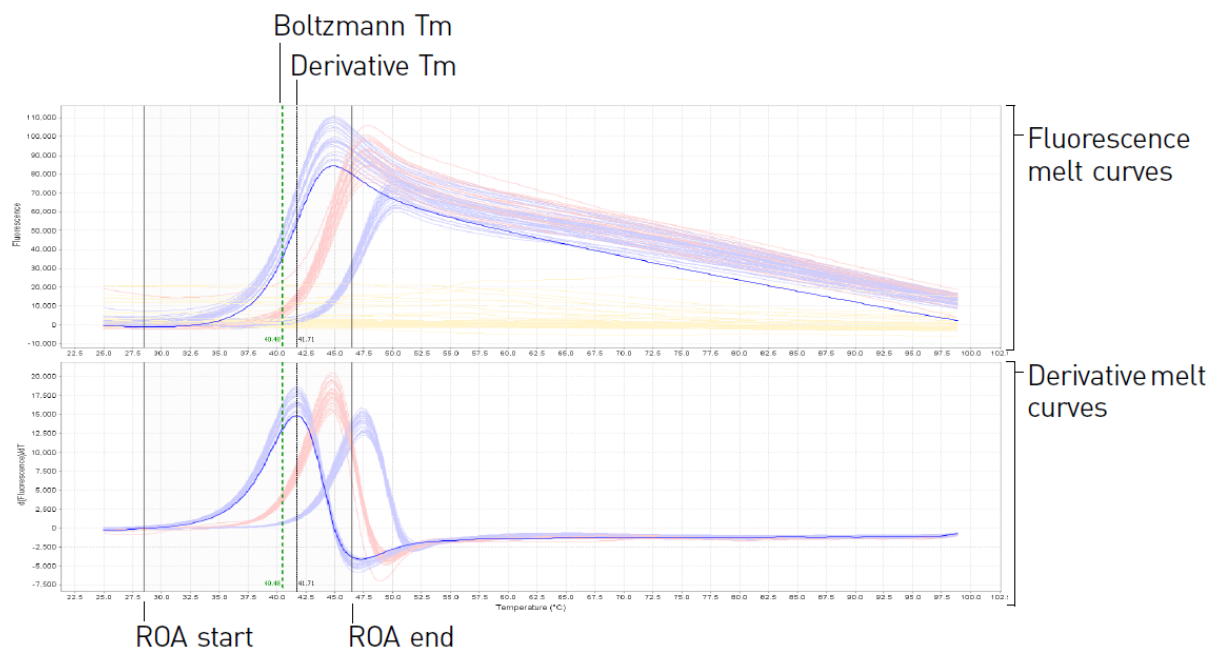
Using the Protein Thermal Shift™ Software, review the melt curves and well table and optimize the analysis in the Well Results screen.

The Well Results screen displays fluorescence and derivative melt curve plots, calculated T_m values, individual well results, and flags. As necessary, edit the analysis settings, edit the baseline, edit the region of analysis, edit the analysis mode, and omit outliers.

This section provides guidance on how to review and interpret the well results. Some common troubleshooting causes are provided. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

About the melt curve plots

In the Well Results screen, the fluorescence data are plotted as fluorescence melt curves and as derivative melt curves. The regions of analysis (ROAs), Boltzmann T_m values, and derivative T_m values are displayed in the melt curve plots.



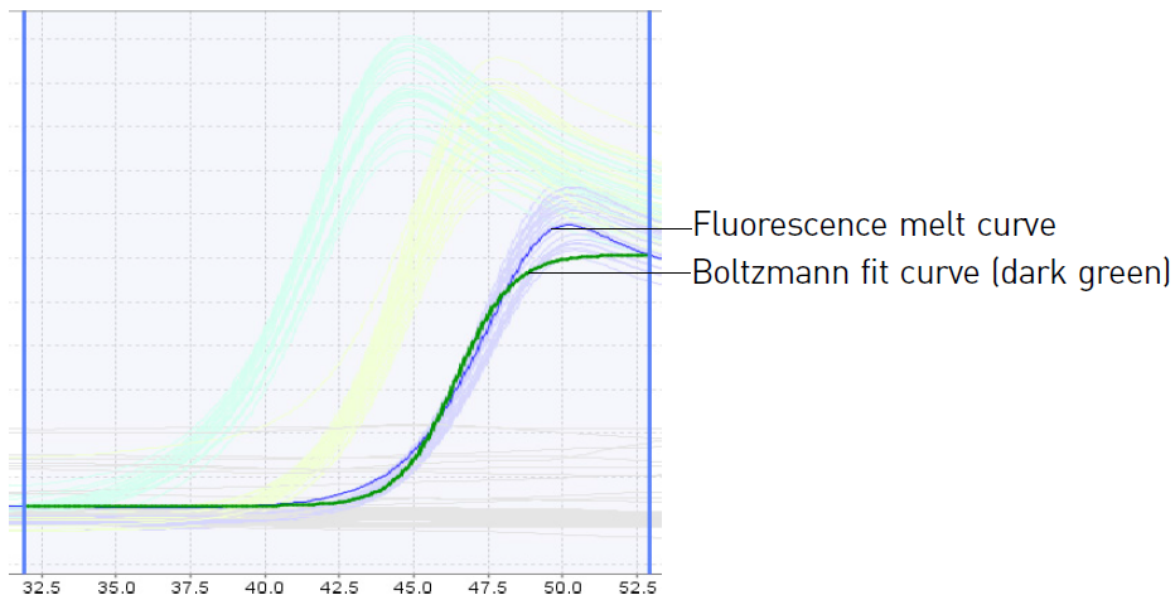
The fluorescence melt curves show noise-reduced fluorescence data from the protein melt curve run. The fluorescence melt curves and the Boltzmann method are used to determine the Boltzmann T_m .

The derivative melt curves are calculated using a first derivative of the fluorescence data at each temperature. The derivative melt curve is used to identify melt phases, or regions of analysis (ROAs), from which the derivative T_m is calculated. In Protein Thermal Shift™ Software, up to six ROAs and derivative T_m values can be automatically detected or manually defined for each melt curve.

Because the Boltzmann T_m and derivative T_m values are calculated independently of each other, the values may be dissimilar.

About the Boltzmann fit curve

The Boltzmann fit curve is calculated for the region of analysis (ROA) according to the Boltzmann equation and is plotted in the fluorescence melt curve plot as a dark green curve.



The Boltzmann fit value corresponds to how similar the fluorescence melt curve and the Boltzmann fit curve are.

Note: When you define the ROA manually, you may observe a gap between the ROA start or end temperature and the start or end of the Boltzmann fit curve. The gap occurs if the defined ROA start or end temperature does not correspond exactly with a fluorescence datapoint.

Review the melt curves

Review the melt curves to visualize the fluorescence and derivative fluorescence data in the example ligand titration study.

Note: For NPC wells, the derivative melt curves are not displayed.

1. In the **Analysis ▶ Well Results** screen, select **Show in Plot**, then select to show these plot components:
 - Unselected Wells
 - Legend
2. Select **Color By**, then select **Ligand** to color the melt curves according to the ligand condition value assigned for the well.

Note: To set the color for each ligand to make the plots easier to distinguish, go to the **Conditions** screen.

3. For each replicate group, select all the wells in the replicate group, then review the fluorescence levels in the melt curves:
- For the NPC wells, do you observe a rise in fluorescence? If so, the wells or protein melt reactions may be contaminated with protein or the dye may interact with a buffer component.
 - For the LOC wells, do you observe a rise in fluorescence? A rise in fluorescence in LOC wells but not in NPC wells may indicate protein contamination in the ligand or ligand-dye interactions.
 - For sample or reference wells, do you observe flat melt curves? If so, condition assignments may be incorrect or a component is missing from the protein melt reactions.
 - Within each replicate group, are the fluorescence melt curves similar to each other? Within each replicate group, are the derivative melt curves similar to each other? If the melt curves for the replicates are dissimilar, pipetting errors may have occurred during reaction setup or condition assignments may be incorrect.

About the analysis mode

In the Protein Thermal Shift™ Software, the analysis mode is the method for defining the regions of analysis (ROA) and determining the derivative T_m for a melt curve.



The analysis mode is displayed in the Analysis Mode column of the Well Table in the Analysis screens.

Analysis mode	Description
Auto: Single T _m	The software detects one melt phase, defines one ROA, and determines one derivative T _m within the ROA. By default, the melt curves are analyzed using the Auto: Single T _m analysis mode.
Auto: Multiple T _m	The software detects more than one melt phase, defines an ROA for each melt phase, and determines one derivative T _m for each ROA. The software can detect up to six ROAs for each melt curve. If the derivative melt curve shows multiple melt phases, change the analysis mode to Auto: Multiple T _m .
Manual	You define the ROA within the software, then the software determines one derivative T _m for each ROA that you defined. You can define up to six ROAs for each melt curve. If you edit or delete an ROA detected by the software, the analysis mode is changed to Manual.

Set the analysis mode

In the example ligand titration study, the melt curves should show one melt phase. If you see more than one peak, the wells or protein melt reactions may be contaminated with protein.


In your own studies, you may observe multiple melt phases. If the derivative melt curves for the replicate group show multiple melt phases, set the analysis mode to Auto: Multiple T_m, then review the derivative melt curves.

1. In the Well Table or in the melt curve plot, select the wells with multiple melt phases.
2. Click  **Auto Analysis Options**, then select **Auto: Multiple T_m**.
3. Click  **Analyze** to reanalyze using the Auto: Multiple T_m analysis mode.
4. Review the number of melt phases (peaks) in the derivative melt curves:
 - Do all replicate groups have the same number of melt phases?
 - For each replicate group, are there outliers with a different number of melt phases than the other samples in the replicate group? You may consider omitting outliers from analysis.

Review the regions of analysis (ROA)

Review the ROAs detected by the software. If necessary, edit the ROAs.

Note: If no melt phases are detected, no ROAs are defined. Negative controls should have no melt phases and no ROAs.

1. In the Analysis > Well Results screen, confirm that each ROA meets the following criteria:
 - For melt curves with one melt phase, the curve within the ROA resembles a sigmoidal profile.
 - At the start temperature, the signal is relatively flat.
 - At the end temperature, the signal has already reached its maximum.
2. For each replicate group, edit the ROAs so that all of the wells in the replicate group have the same ROA:
 - a. Select the replicates, click  (Define ROA) in the toolbar above the melt curve plots, then click-drag an area in one of the plots to define a melt phase and replace the ROA. Repeat for each melt phase you identify.
 - b. If necessary, adjust the start and end temperatures for each ROA:
 - **To move the ROA:** Starting from within the ROA, click-drag the ROA.
 - **To adjust the start and end temperatures individually:** Click-drag the ROA start or the end line.



IMPORTANT! Make sure that the fluorescence at the start temperature is lower than the fluorescence at the end temperature.

3. Click  **Analyze** to reanalyze using the edited ROAs.

Note: After you edit ROAs, the analysis mode changes from Auto to Manual.

Review the flags and Tm values


Review the flags and Tm values, then consider editing the analysis settings and/or omitting wells before you review the replicate results.

1. In the **Well Results** screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann Tm
 - Derivative Tm
 - Unselected Wells
 - Legend
2. Review flagged wells in the Well Table:
 - a. Click the  (Flag Indicator) column header to sort the wells according to the number of flags applied to the well.
 - b. Scroll the table to the right to view the flags that are applied to the wells.




- c. For flagged wells, select the well in the Well Table, then review the melt curves for the well, compared to the other wells in the replicate group.
 - d. Omit wells from analysis as necessary.
3. For each replicate group, select the wells in the replicate group, then review the Tm B (Boltzmann Tm) and the Tm D (derivative Tm) in the Well Table and in the melt curves:

Note: In the melt curves, the Boltzmann Tm is a green dashed vertical line, and the derivative Tm is black dotted vertical line.

- Are the Tm B or Tm D values significantly different from the Tm values for other wells in the replicate group? Do any replicates have melt curves that are inconsistent with the other melt curves for the replicate group? If so, you may consider omitting wells from analysis.
 - Are the melt curves within the replicate group similar?
4. If you omitted any wells from the analysis, click  **Analyze**.

Review the Boltzmann fit

1. In the **Well Results** screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann fit
 - Unselected Wells
 - Legend

2. Scroll through each well in the Well Table, then compare the fluorescence melt curve to the Boltzmann fit curve (dark green thick curve) and review the value in the B Fit (Boltzmann fit) column of the Well Table:
 - How well does the fluorescence melt curve correspond to the Boltzmann fit curve?
 - Is the B Fit value close to 1?
 - Is the B Fit value similar among wells in the replicate group?

Note: When you define the ROA manually, you may observe a gap between the ROA start or end temperature and the start or end of the Boltzmann fit curve. The gap occurs if the defined ROA start or end temperature does not correspond exactly with a fluorescence datapoint.

Example well results

For the ViiA™ 7 System example ligand titration study using the Protein Thermal Shift™ Control Protein and the Protein Thermal Shift™ Control Ligand, observe the following:

- Fluorescence levels are flat in the NPC and LOC wells.
- For the sample and reference replicate groups, the fluorescence levels as displayed in the fluorescence melt curve are not significantly different, so you do not need to edit the baseline.
- The sample and reference wells contain one peak in the derivative melt curve.
- For each replicate group, the Boltzmann T_m values are similar and there are no outliers.
- For each replicate group, the derivative T_m values are similar and there are no outliers.
- For each well, the region of analysis meets the recommended criteria.

Review the replicate results

Using the Protein Thermal Shift™ Software, review the T_m statistics for the replicate groups and look for positive hits in the **Analysis > Replicate Results** screen.

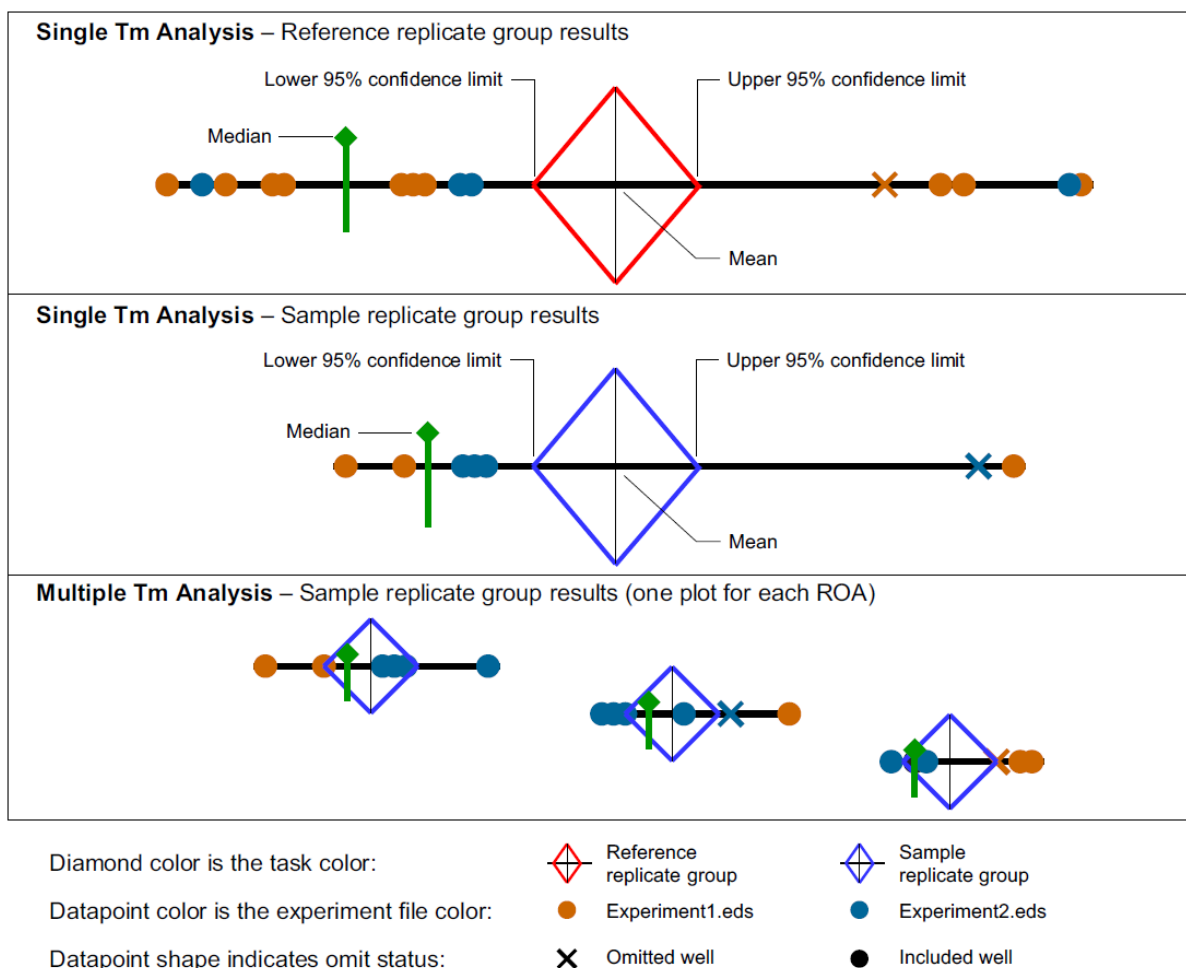
This section provides guidance on how to review and interpret the replicate results. Some common troubleshooting causes are provided. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

About the T_m statistics

The ΔT_m values are calculated for each ROA by subtracting the mean T_m for the replicate group from the mean T_m for the reference replicate group. If there is more than one ROA, the ΔT_m values are calculated for each ROA in the order that they appear (left to right).

About the replicate results plot


The Replicate Results Plot shows T_m statistics for each replicate group and T_m datapoints for each well in the replicate group. You can select to view the Boltzmann T_m statistics, derivative T_m statistics, Δ Boltzmann T_m statistics, or Δ Derivative T_m statistics. The plot shows the calculated T_m or ΔT_m value for each sample in the replicate group. The mean T_m or ΔT_m value and the upper and lower 95% confidence limits are identified by colored diamonds - red diamonds for the reference replicate groups and blue diamonds for the sample replicate groups.



Review the boltzmann T_m statistics

In the **Analysis ▶ Replicate Results** screen, review the Boltzmann T_m statistics to evaluate the variability among replicates. The Boltzmann T_m statistics are plotted along the x-axis for each replicate group.


For the example ligand titration study, we provide instructions for reviewing the Boltzmann T_m statistics. For your own studies, you may also review derivative T_m statistics.

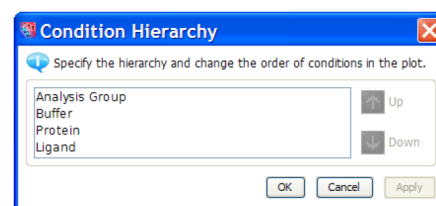
1. In the **Analysis ▶ Replicate Results** screen, select  **Plot by**, then select **T_m-Boltzmann**.

Note: Select **T_m-Derivative** to review the derivative T_m statistics.



2. Specify the condition hierarchy to group the replicate plots and change the order of conditions in the plot:

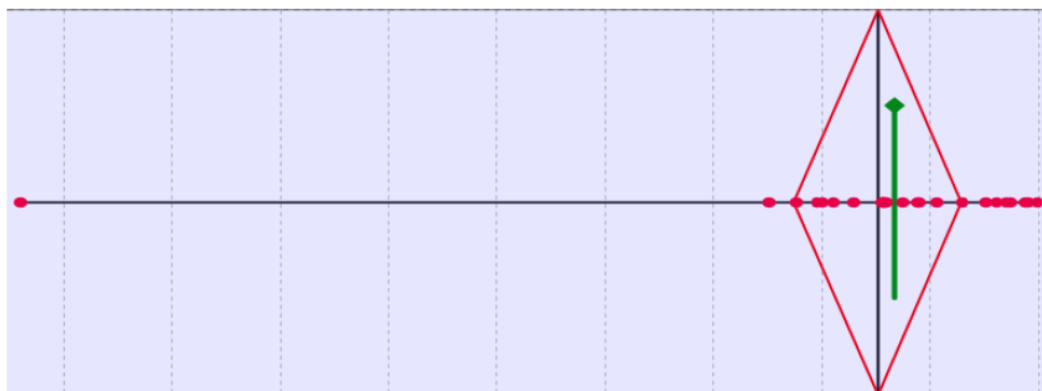
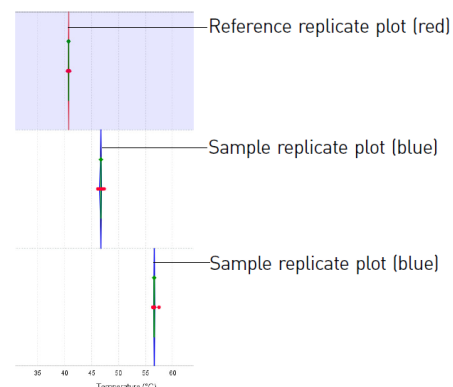
Note: Changing the condition hierarchy does not affect the results; it only changes how the replicate plots are grouped and the order in which the conditions are displayed in the Replicate Results Plot.

- a. Click  **Condition Hierarchy** above the top right corner of the plot.
- b. In the dialog box, select a condition, then use the Up and Down arrows to change the hierarchy of conditions.
- c. For the example ligand titration study, continue to move the conditions up and down until you obtain this hierarchy:
The condition at the top-most level of the hierarchy is displayed on the far-right side of the Replicate Results Plot:

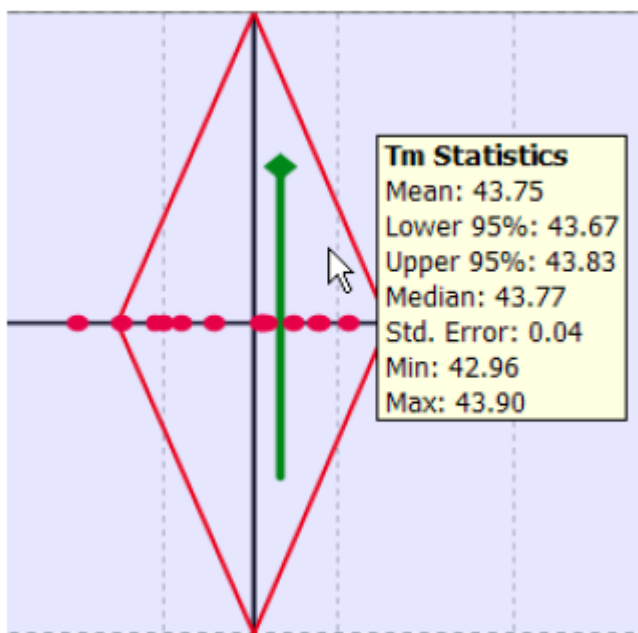


Ligand Protein Buffer Analysis Group

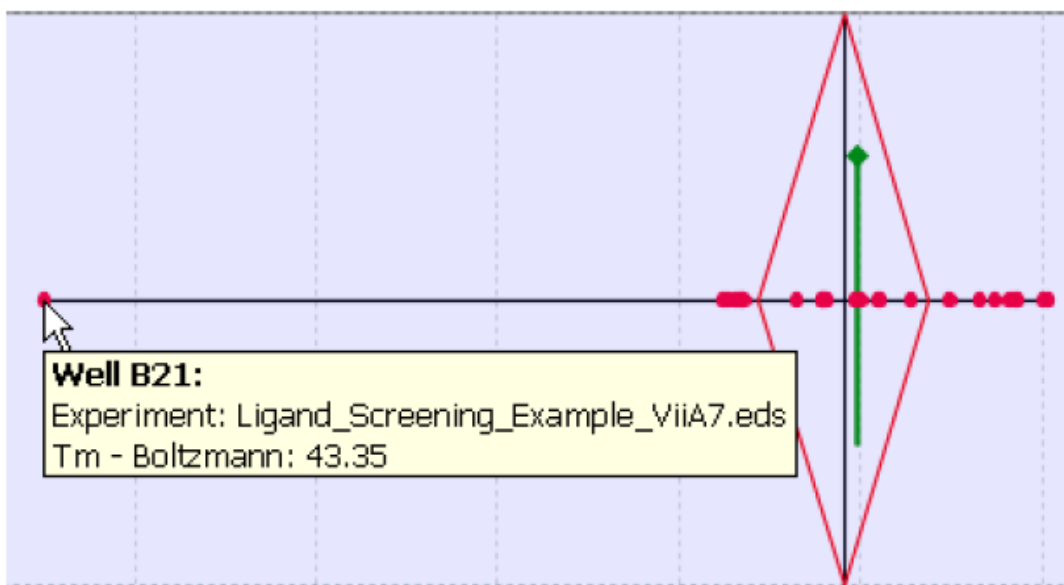
3. Scan the Replicate Results Plot to review the conditions that affect the T_m values relative to the reference replicate group.
4. Review the T_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot one or more times.
 - b. Move the plot: Click , then click-drag the plot until the replicate plot of interest is in view.




- c. Place the cursor within the diamond, then wait to view a tooltip with the T_m statistics for the replicate group.



- d. To examine outliers, place the cursor over a datapoint, then wait to view a tooltip with the well information, experiment file name, and the T_m selected for the plot.



Note: Click  to restore the default zoom.

5. In the Replicate Groups table, review the Boltzmann T_m statistics for each replicate group:
 - **T_m B - Std. Error** (standard error of the mean for the Boltzmann T_m): Is the value low? If the value is high, review the data for each replicate.
 - **T_m B - Min** and **T_m B - Max™** (minimum and maximum Boltzmann T_m values for the replicate group): Is the range of Boltzmann T_m values for the replicate group within 1 degree? If the range of T_m values exceeds 1 degree, review the data for each replicate.
6. Omit outliers as necessary, then click  **Analyze**.



Review the positive hits

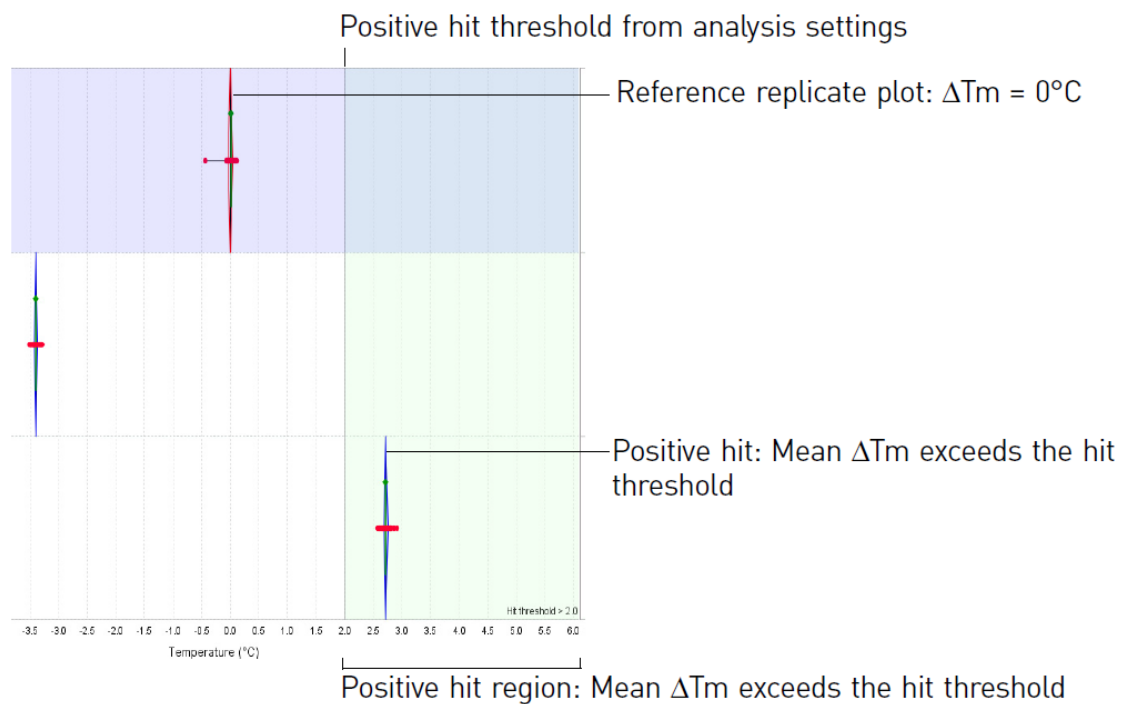
In the Analysis > Replicate Results screen, review the positive hits to identify the conditions that produce the maximum effect on thermal stability. Replicate groups with positive hits have ΔT_m values that exceed the threshold set in the analysis settings.



Note: The positive hits are determined according to the analysis settings that you specified (“Review the analysis settings” on page 42).

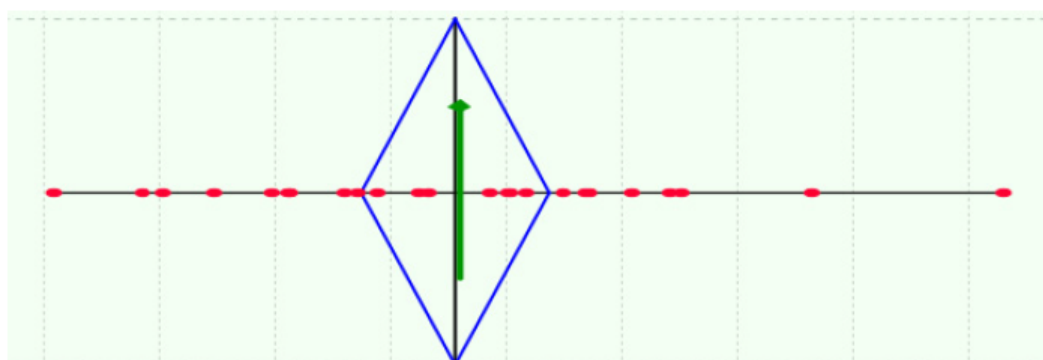
For the example ligand titration study, review the ΔT_m -Boltzmann positive hits. In your own studies, you may choose to review the ΔT_m - Derivative positive hits.

Note: You must specify the reference replicate group to calculate ΔT_m values and to determine positive hits.

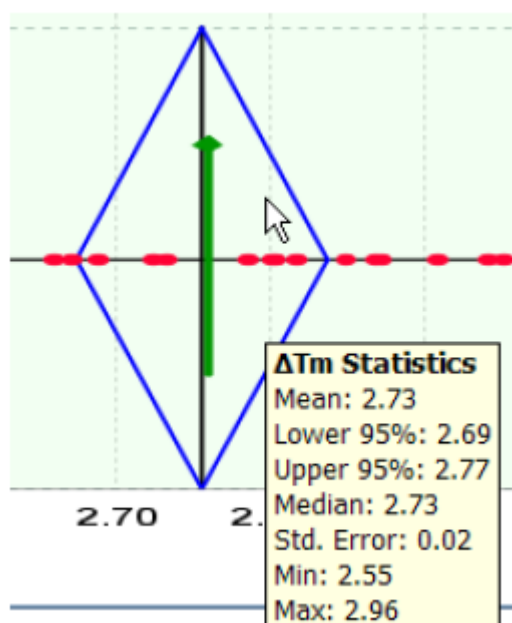
1. In the Analysis > Replicate Results screen, select the  **Plot by** menu, then select **ΔT_m - Boltzmann**.
2. Select the  **Show in Plot** menu, then select to show **Positive Hits**.
3. Scan the Replicate Results Plot for positive hits.
The positive hits are replicate plot(s) within the green shaded area and labeled with the hit threshold in the bottom right corner of the plot.



4. Review the ΔT_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot as many times as you need to magnify the plot.
 - b. Move the plot: Click , then click-drag the plot until the replicate plot you want to review is in view.



- c. Place the cursor over the replicate plot, then wait to view a tooltip with the ΔT_m statistics for the replicate group.



Note: Click to restore the default zoom.

5. In the Replicate Groups table, review the positive hits (🟢) in the Hits B column and review the ΔT_m B statistics in the table.

Example replicate results

For the ViiA™ 7 System example ligand titration study using the starter kit, observe the following:

- The control ligand stabilizes the protein in a concentration-dependent manner.
- The protein melt reactions that contained 1.0 mM control ligand increased the thermal stability of the control protein beyond the threshold set in the analysis settings (ΔT_m Boltzmann > 2.0°C).

Next steps

The use of Protein Thermal Shift™ studies to determine the relative thermal stability of proteins (ΔT_m values) is precise and consistent with analysis using calorimetry or spectroscopy methods. After you perform Protein Thermal Shift™ studies to screen for and identify positive hits, you can validate the positive hits using calorimetry or spectroscopy methods to determine the thermodynamic stability.

5

Buffer screening studies

Perform a buffer screening study to identify the optimal buffer, pH, or salt concentration for storing a protein. Buffer screening may also be used to identify a buffer system for a protein that aggregates or precipitates from solution or to improve protein crystal formation for x-ray crystallography.

“Create and set up an experiment file for the instrument run” on page 56



“Prepare the protein melt reactions” on page 59



“Run the protein melt reactions” on page 60



“Set up the analysis” on page 62



“Review the well results” on page 66



“Review the replicate results” on page 71

Example experiment files and plate template files are located in the examples folder: `...Program Files\Applied Biosystems\Protein Thermal Shift Software\examples`, where `...` is the installation drive.

To view the data for the buffer screening example used in this chapter, use the `Buffer_Screening_Example_ViiA7.eds` and `Buffer_Screening_Example_Setup_ViiA7.csv` files in the `ViiA7 Example Files` folder.

General guidelines

For general guidelines that apply to all Protein Thermal Shift™ studies, see “General guidelines” on page 12.

Experimental design

With Protein Thermal Shift™ studies, you can perform a series of buffer screening studies to identify the optimal buffer, pH, and salt concentration for a protein. You may also perform a buffer screening study before performing a mutation screening study or a ligand screening study.

Replicates and controls

For buffer screening studies, we recommend that you prepare:

- At least 4 replicates of each reaction
- At least 4 replicates of no protein controls (NPCs)

Create and set up an experiment file for the instrument run

This section contains general settings for creating and setting up an experiment file. For detailed instructions, see the following sections.

Instrument	Page
QuantStudio™ 7 Pro Real-Time PCR System QuantStudio™ 6 Pro Real-Time PCR Instrument	page 15
QuantStudio™ 5 Real-Time PCR System QuantStudio™ 3 Real-Time PCR System QuantStudio™ 1 Real-Time PCR System	page 18
QuantStudio™ 7 Flex Real-Time PCR System QuantStudio™ 6 Flex Real-Time PCR System QuantStudio™ 12K Flex Real-Time PCR System	page 22
ViiA™ 7 Real-Time PCR System	page 28
StepOne™ Real-Time PCR System StepOnePlus™ Real-Time PCR System	page 31
7500 Real-Time PCR System 7500 Fast Real-Time PCR System	page 34

Experiment file settings

Setup	Setting
Experiment properties	<ul style="list-style-type: none"> • Experiment type: Melt Curve • Reagents: Other • Ramp speed: Fast or Standard
Target properties	<ul style="list-style-type: none"> • Reporter: ROX™ • Quencher: None
Plate layout	<ul style="list-style-type: none"> • Assign targets to all wells in use • Passive reference: None

(continued)

Setup	Setting
Run method	<ul style="list-style-type: none"> • Reaction Volume Per Well: 20 μL • Thermal profile: <ul style="list-style-type: none"> – Step 1, Temp: 25°C, Time: 2 minutes – Step 2, Temp: 99°C, Time: 2 minutes • Ramp mode: Continuous • Ramp rate: Step 1: 1.6°C/s, Step 2: 0.05°C/s for the following systems. <ul style="list-style-type: none"> – QuantStudio™ 7 Pro Real-Time PCR System – QuantStudio™ 6 Pro Real-Time PCR System – QuantStudio™ 5 Real-Time PCR System – QuantStudio™ 3 Real-Time PCR System – QuantStudio™ 1 Real-Time PCR System – QuantStudio™ 7 Flex Real-Time PCR System – QuantStudio™ 6 Flex Real-Time PCR System – QuantStudio™ 12K Flex Real-Time PCR System – ViiA™ 7 Real-Time PCR System Ramp rate: 1% for the following systems. <ul style="list-style-type: none"> – StepOne™ Real-Time PCR System – StepOnePlus™ Real-Time PCR System – 7500 Real-Time PCR System – 7500 Fast Real-Time PCR System • Optical Filters (ViiA™ 7 System only): <ul style="list-style-type: none"> – Coupled filters: x2(520\pm10)—m2(558\pm11) – Decoupled filters: x1(470\pm15)—m3(586\pm10)

Defining and assigning targets

We recommend that you define and assign targets so you can review the melt curves for the replicate groups in the Real-Time PCR System Software before importing the experiment files into the Protein Thermal Shift™ study. Define the following targets:

- One target for each buffer condition
- Define a target for the no protein control (NPC)

IMPORTANT! Make sure that the well assignments correspond exactly with the arrangement of reactions in the reaction plate. Setup errors may result in an incorrect grouping of replicates.




















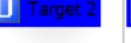

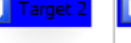





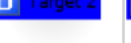




































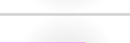
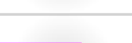

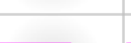
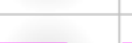











Example targets

In the ViiA™ 7 Real-Time PCR System buffer screening example file, targets were defined for each buffer and the no protein control (NPC).

Target Name	Reporter	Quencher	Color
Target 1	ROX	None	Red
Target 2	ROX	None	Blue
Target 3	ROX	None	Green
Target 4	ROX	None	Cyan
NPC	ROX	None	Magenta

Example plate layout

In this example, columns 9–24 and rows K–P are not in use.

	1	2	3	4	5	6	7	8
A	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1
B	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1	 Target 1
C	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2
D	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2	 Target 2
E	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3
F	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3	 Target 3
G	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4
H	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4	 Target 4
I	 NPC	 NPC	 NPC	 NPC	 NPC	 NPC	 NPC	 NPC
J	 NPC	 NPC	 NPC	 NPC	 NPC	 NPC	 NPC	 NPC

Prepare the protein melt reactions

Protein melt reaction stability

For consistency in T_m values, we recommend that you keep the protein melt reactions on ice until you are ready to load the instrument and start the run.

If the protein is thermally stable at ambient temperatures, you may consider preparing the reactions in advance and leave the reaction plate at ambient temperature, protected from light. However, the fluorescence levels will decrease over time and the T_m values will vary, depending on the protein and its thermal stability. If you want to prepare the reaction plates in advance, we recommend that you first determine the benchtop stability of your protein melt reactions.

Required materials

Required materials for protein melt reactions:

- Protein Thermal Shift™ Dye (1000X)
- Protein Thermal Shift™ Buffer
- Water
- Protein
- Buffers and/or buffer components
- MicroAmp™ Optical Reaction Plate appropriate for your real-time PCR instrument
- MicroAmp™ Optical Adhesive Film appropriate for your reaction plate

Prepare the protein melt reactions

We recommend that you prepare four replicates of each reaction.

1. Prepare a fresh dilution of Protein Thermal Shift™ Dye (1000X) to 8X.
2. Place the appropriate reaction plate or tubes on ice, then prepare the protein melt reactions:
 - Make sure that the arrangement of reactions in the reaction plate corresponds exactly with the well assignments in the experiment file.

IMPORTANT! Setup errors may result in an incorrect grouping of the data and incorrect T_m statistics.

- Add reaction components to the plate in the order listed.

Component	Volume
Protein Thermal Shift™ Buffer	5.0 µL
Water + protein + buffer and/or buffer components	12.5 µL
Diluted Protein Thermal Shift™ Dye (8X)	2.5 µL
Total volume for each control reaction	20.0 µL

3. Pipet each reaction up and down 10 times to mix well.
4. Seal the plate with MicroAmp™ Optical Adhesive Film, centrifuge at 1000 rpm for 1 minute, then place it on ice.

Run the protein melt reactions

Load and run the protein melt reactions on a supported Applied Biosystems™ Real-Time PCR System, then analyze and save the experiment file before importing the experiment file into the Protein Thermal Shift™ Software.

IMPORTANT! Keep the protein melt reactions on ice until you load the instrument.

Load and run the reactions

1. Transfer the experiment file that you created for the run to the computer that is connected to the instrument, if needed.
2. In the home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file you created for the run.
3. Load the reaction plate into the instrument.
4. In the software, start the run.

Review the melt curves

Using the Real-Time PCR System Software, open the experiment file from the completed instrument run, analyze and save the experiment file, then review the melt curves.

Note: You must analyze and save the experiment file in the Real-Time PCR System Software before you can import it into the Protein Thermal Shift™ Software.

Some common troubleshooting causes are provided here. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

1. In the Home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file from the instrument run.
2. View the melt curves:

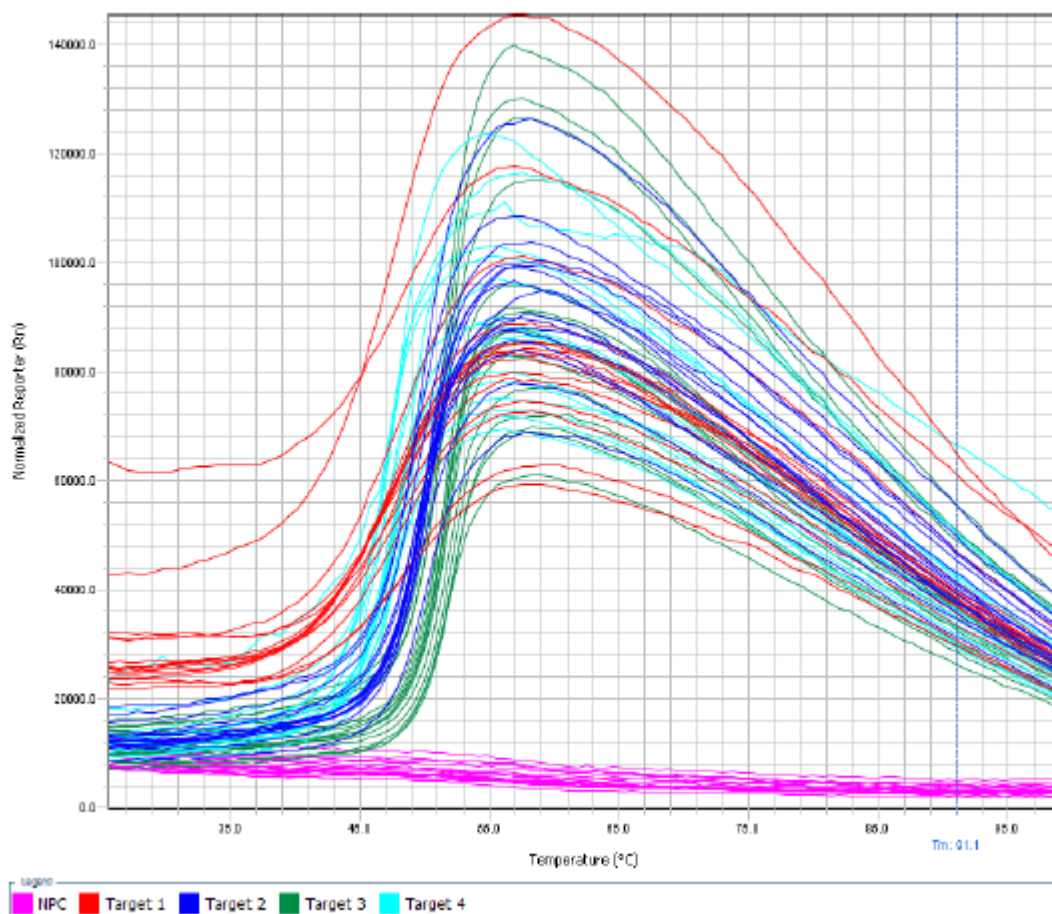
Real-time PCR System Software	View the melt curve
ViiA™ 7 Software	Click Analysis ▶ Melt Curve Plot in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.
StepOne™ Software or 7500 Software	Click Analysis ▶ Melt Curve in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.

3. Review the melt curves:
 - Do you see fluorescence signals in all of the sample wells?
No fluorescence signals in the sample wells may indicate missing dye or protein or an instrument problem.
 - Do you see flat fluorescence levels in the NPC wells?
High fluorescence levels in the NPC wells may indicate protein contamination in the wells or protein melt reactions; or it may indicate that the dye interacts with a component in the buffer.
 - Do the replicates have similar melt curves?
4. Save, then close the experiment file.

Note: The melt curves in the real-time PCR software may not exactly match the melt curves in the Protein Thermal Shift™ Software. When the experiment files are imported into the Protein Thermal Shift™ Software, the Protein Thermal Shift™ Software reduces the noise in the fluorescence data.

Example melt curves

Melt curves for the buffer screening example file from the ViiA™ 7 Real-Time PCR System:





Set up the analysis

This section provides instructions for setting up the Protein Thermal Shift™ study using Protein Thermal Shift™ Software v1.4.

Setup guidelines

- The experiment files that you import into the study must contain analyzed melt curve data from a complete melt curve run.
- Set up the analysis group so that it contains experiment files from only one instrument.

Create and set up the study

1. In the Home screen of the Protein Thermal Shift™ Software, click  **Create Study**.
2. Complete the **Setup > Properties** screen:
 - The Study Name cannot be more than 100 characters and cannot contain these characters: / \ * " ? < > | . ,
 - The instrument selection must match the instrument type that you used to run the protein melt reactions and generate the experiment files.
3. In the **Setup > Conditions** screen, define the conditions and condition values, then list the analysis groups for your study.
 - a. In the **Condition Names** pane, click  **Add** to add up to 20 custom parameters for the study, such as pH.
 - b. Select a condition of interest, then in the pane for the values for the condition, add values for the condition.













Condition Names		Values for Salt					
 Add		 Add		 Import	 Up		Down
Name		Name	Color	Description	Comments	# of Wells Assigned	
Protein		0 mM					0
Ligand		50 mM					0
Buffer		100 mM					0
Salt		150 mM					0
pH							

Figure 1 Example with 4 concentrations for Salt added.

- c. In the **Analysis Groups** pane, create groups, then assign wells to each group to perform a more focused analysis with the software, if needed.
4. In the **Setup > Experiment Files** screen, click  **Import**, then select the experiment file (EDS) for the instrument type that you selected for the study. Repeat for each experiment file to import into the study.

Note: The experiment file name and data are imported. Plate setup information is not imported.



5. For each experiment file, assign the conditions, task, and analysis group to the wells that contained protein melt reactions.
 - a. In the **Setup > Experiment Files** screen, click-drag wells in the plate to select multiple wells. If the wells are not continuous, ctrl+click the wells of interest.
 - b. Right-click one of the selected wells, then click **Assign Well Content**.
 - c. In the dialog box that appears, select the value for each condition that applies to the selected wells, then click **Done**.
In the **Task** dropdown list, you can select **Sample**, **Reference** or **NPC** to assign the well as an unknown sample, a reference sample, or a no protein control.

- d. Repeat step 5 and substep 5a for each condition.

You can click **Auto-Fill Settings** to assign the wells faster.

6. Ensure that the condition assignments correspond exactly with the contents of the reaction plate.

IMPORTANT! Setup errors can result in an incorrect grouping of the data and incorrect T_m statistics.

7. For each analysis group, assign the reference replicate group:
 - a. Select the wells for the replicate group to use as the reference.
 - b. Click  **Assign**, then select **Reference** as the Task.
8. Click  **Save** in the toolbar to save and analyze the study.

Assign wells automatically

You can use auto fill settings to assign wells faster.

Alternatively, you can click **Load Plate Template** to import a Microsoft™ Excel™ spreadsheet with the well information.

Create and set up the study.

1. In the **Setup > Experiment Files** screen, assign one or more conditions to one or more rows or columns of wells.
2. Click **Auto-Fill Settings ▶ Dock**.
3. In the **Auto-Fill Settings** pane, make the following selections for each condition.
 - a. Click the ✓ (**checkmark**) to deselect a condition so that values are not automatically assigned for the condition, if needed.
 - b. For each selected condition, select **Copy** to copy the value for the condition, or select **Series** to extend the sequence for the series of values for the condition.



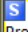
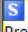
For each series, the values are added in the order in which the values were created in the study setup.
4. For each column or row that you want to fill, click-drag to select wells from which the condition values are copied or extended.

5. Place the pointer over the last selected well of the column or row, then drag the pointer over the wells that you want to fill.

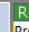
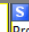
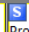
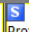
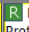
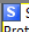

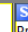
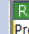
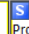
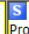
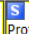
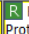
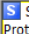
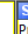
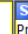
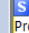
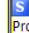

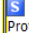
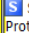
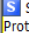
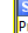
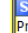
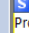
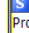
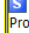
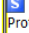
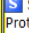
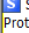
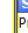
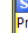
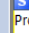
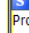
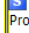
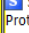
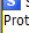
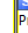
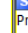
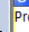
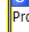
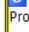


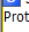
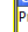

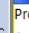
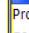
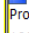
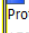
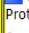
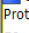
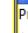
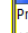
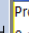
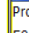
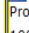
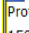
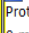
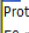
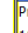
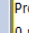
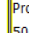
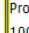
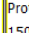
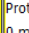
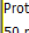
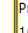
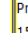
6. Click  **Save** to save the study.

The remaining wells are filled according to the selections.

Example plate setup

7	8	9	10
 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1		
 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1		


Well 18
Task: Sample
Analysis Group: AG 1
Protein: Protein A
Salt: 150 mM

	1	2	3	4	5	6	7	8
A	 Reference Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Reference Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
B	 Reference Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Reference Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
C	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
D	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
E	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
F	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
G	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
H	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1
I	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1	 Sample Protein A 0 mM AG 1	 Sample Protein A 50 mM AG 1	 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1

Review the analysis settings

Review and edit the analysis settings to optimize the analysis for your study.

For the examples shown in this user guide, no changes were made to the analysis settings. If you are reviewing the data in the example studies provided with the software, try revising the analysis settings to see how the settings affect the positive hits and the flags.

1. Click  **Analysis Settings** in the toolbar.
2. On the Positive Hit tab, specify the ΔT_m -Boltzmann and ΔT_m -Derivative values to indicate a positive hit:
 - Select **>** to identify buffer conditions that increase protein thermal stability, or select **<** to identify buffer conditions that decrease protein thermal stability.
 - Enter the number of degrees ($^{\circ}\text{C}$) of T_m shift relative to the reference to indicate a positive hit.
3. On the Flags tab, specify settings for applying flags:
 - a. Select the flags to use in the analysis.
 - b. For the High Background, High NPC, Low Signal, and Poor Fit flags, specify the condition and threshold for applying the flag.
4. Apply the analysis settings and analyze:
 - Click **Apply** to apply the analysis settings and reanalyze while keeping the Analysis Settings dialog box open.
or
 - Click **OK** to apply the analysis settings, reanalyze, and close the Analysis Settings.

Review the well results

Using the Protein Thermal Shift™ Software, review the melt curves and well table and optimize the analysis in the Analysis > Well Results screen.



The Well Results screen displays fluorescence and derivative melt curve plots, calculated T_m values, individual well results, and flags. As necessary, edit the analysis settings, edit the baseline, edit the region of analysis, edit the analysis mode, and omit outliers.

This section provides guidance on how to review and interpret the well results. Some common troubleshooting causes are provided. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

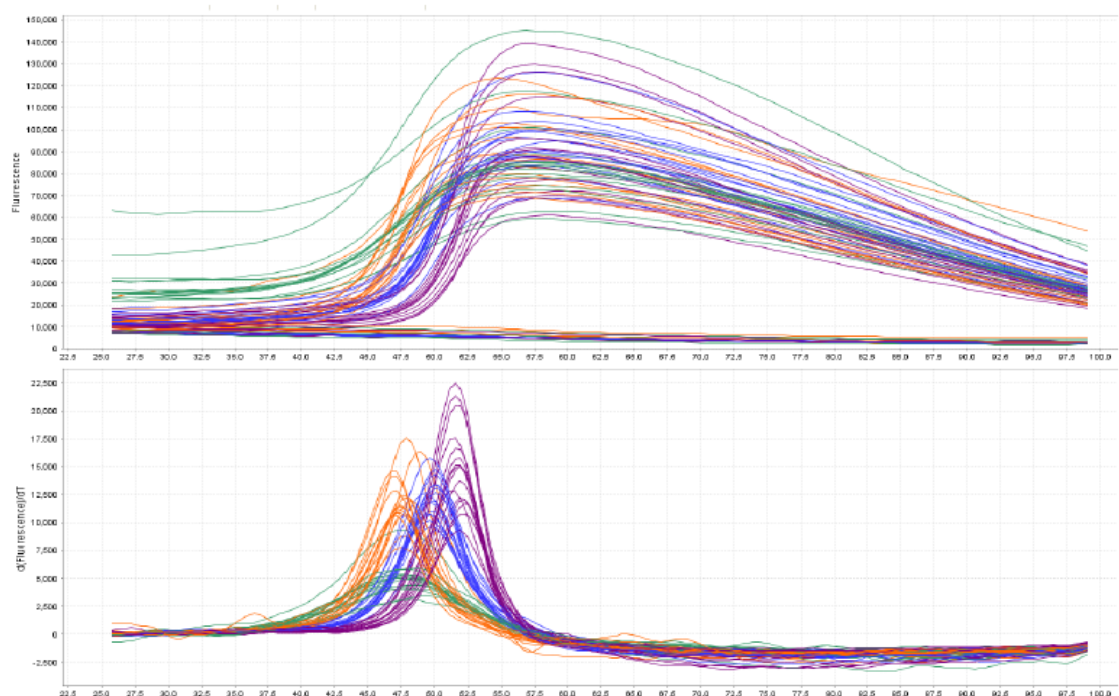
Review the melt curves



Review the melt curve plots to visualize the fluorescence and derivative fluorescence data. If necessary, change the analysis mode.

Note: For NPC wells, the derivative melt curves are not displayed.

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Unselected Wells
 - Legend
2. select  **Color By**, then select a buffer condition (for example, **Buffer** or **Salt**) to color the melt curves according to the buffer condition value assigned for the well.

Note: To set the color for each buffer condition to make the plots easier to distinguish, go to the Setup > Conditions screen.




3. For each replicate group, select all the wells in the replicate group, then review the fluorescence levels in the melt curves:
 - For the NPC wells, do you observe a rise in fluorescence? If so, the wells or protein melt reactions may be contaminated with protein or the dye may interact with a buffer component.
 - For sample or reference wells, do you observe flat melt curves? If so, condition assignments may be incorrect or a component is missing from the protein melt reactions.
 - Within each replicate group, are the fluorescence melt curves similar to each other? Within each replicate group, are the derivative melt curves similar to each other? If the melt curves for the replicates are dissimilar, pipetting errors may have occurred during reaction setup or condition assignments may be incorrect.
4. If the derivative melt curves for the replicate group show multiple melt phases, set the analysis mode to Auto: Multiple T_m, then review the derivative melt curves:
 - a. In the Well Table or in the melt curve plot, select the wells with multiple melt phases.
 - b. Click  **Auto Analysis Options**, then select **Auto: Multiple T_m**.
 - c. Click  **Analyze** to reanalyze using the Auto: Multiple T_m analysis mode.
 - d. Review the number of melt phases in the derivative melt curves:
 - Do all replicate groups have the same number of melt phases?
 - For each replicate group, are there outliers with a different number of melt phases than the other samples in the replicate group? You may consider omitting outliers from analysis.

Review the regions of analysis (ROA)

Review the ROAs detected by the software. If necessary, edit the ROAs.

Note: If no melt phases are detected, no ROAs are defined. Negative controls should have no melt phases and no ROAs.

1. In the Analysis > Well Results screen, confirm that each ROA meets the following criteria:
 - For melt curves with one melt phase, the curve within the ROA resembles a sigmoidal profile.
 - At the start temperature, the signal is relatively flat.
 - At the end temperature, the signal has already reached its maximum.
2. For each replicate group, edit the ROAs so that all of the wells in the replicate group have the same ROAs:
 - a. Select the replicates, Click  (Define ROA) in the toolbar above the melt curve plots, then click-drag an area in one of the plots to define a melt phase and replace the ROA. Repeat for each melt phase you identify.

- b. If necessary, adjust the start and end temperatures for each ROA:
 - **To move the ROA:** Starting from within the ROA, click-drag the ROA.
 - **To adjust the start and end temperatures individually:** Click-drag the ROA start or the end line.



IMPORTANT! Make sure that the fluorescence at the start temperature is lower than the fluorescence at the end temperature.

3. Click  **Analyze** to reanalyze using the edited ROAs.

Note: After you edit ROAs, the analysis mode changes from Auto to Manual.

Review the flags and Tm values

Review the flags and Tm values, then consider editing the analysis settings and/or omitting wells before you review the replicate results.

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann Tm
 - Derivative Tm
 - Unselected Wells
 - Legend
2. Review flagged wells in the Well Table:
 - a. Click the  (Flag Indicator) column header to sort the wells according to the number of flags applied to the well.
 - b. Scroll the table to the right to view the flags that are applied to the wells.



- c. For flagged wells, select the well in the Well Table, then review the melt curves for the well, compared to the other wells in the replicate group.
- d. Omit wells from analysis as necessary.


3. For each replicate group, select all the wells in the replicate group, then review the Tm B (Boltzmann Tm) and the Tm D (derivative Tm) in the Well Table and in the melt curves:

Note: In the melt curves, the Boltzmann Tm is a green dashed vertical line, and the derivative Tm is black dotted vertical line.

- Are the Tm B or Tm D values significantly different from the Tm values for other wells in the replicate group? Do any replicates have melt curves that are inconsistent with the other melt curves for the replicate group? If so, you may consider omitting wells from analysis.
- Are the melt curves within the replicate group similar?

4. If you omitted any wells from the analysis, click ► **Analyze**.

Review the Boltzmann fit

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann fit
 - Unselected Wells
 - Legend
2. Scroll through each well in the Well Table, then compare the fluorescence melt curve to the Boltzmann fit curve (dark green thick curve) and review the value in the B Fit (Boltzmann fit) column of the Well Table:
 - How well does the fluorescence melt curve correspond to the Boltzmann fit curve?
 - Is the B Fit value close to 1?
 - Is the B Fit value similar among wells in the replicate group?

Note: When you define the ROA manually, you may observe a gap between the ROA start or end temperature and the start or end of the Boltzmann fit curve. The gap occurs if the defined ROA start or end temperature does not correspond exactly with a fluorescence datapoint.

Example well results

For the ViiA™ 7 Real-Time PCR System example, observe the following:

- Fluorescence levels are flat in the NPC wells.
- For the sample and reference replicate groups, the fluorescence levels as displayed in the fluorescence melt curve are not significantly different, so you do not need to edit the baseline.
- The sample and reference wells contain one peak in the derivative melt curve.
- Within each replicate group:
 - The fluorescence melt curves and the Boltzmann Tm values are similar.
 - The derivative melt curves and the derivative Tm values are similar and there are no outliers.
- For each well, the region of analysis meets the recommended criteria.


Review the replicate results

Using the Protein Thermal Shift™ Software, review the T_m statistics for the replicate groups and look for positive hits in the Analysis > Replicate Results screen.

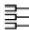
This section provides guidance on how to review and interpret the replicate results. Some common troubleshooting causes are provided. For more detailed information on troubleshooting, see Chapter 8, “Troubleshooting”.

Review the T_m statistics



In the Analysis > Replicate Results screen, review the T_m statistics to evaluate the variability among replicates. The T_m statistics for each ROA are plotted along the x-axis for each replicate group.

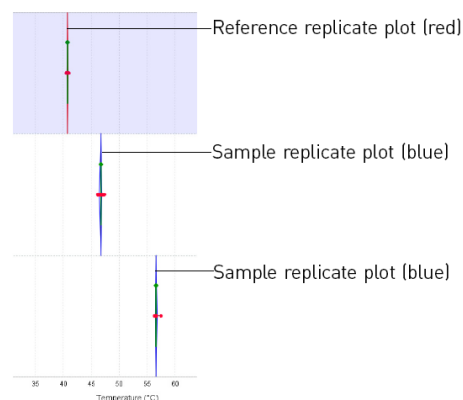
1. In the Analysis > Replicate Results screen, select  **Plot by**, then select the type of T_m statistics to review:
 - **T_m-Boltzmann**
 - **T_m-Derivative**
2. Specify the condition hierarchy to group the replicate plots and change the order of conditions in the plot:

Note: Changing the condition hierarchy does not affect the results; it only changes how the replicate plots are grouped and the order in which the conditions are displayed in the Replicate Results Plot.

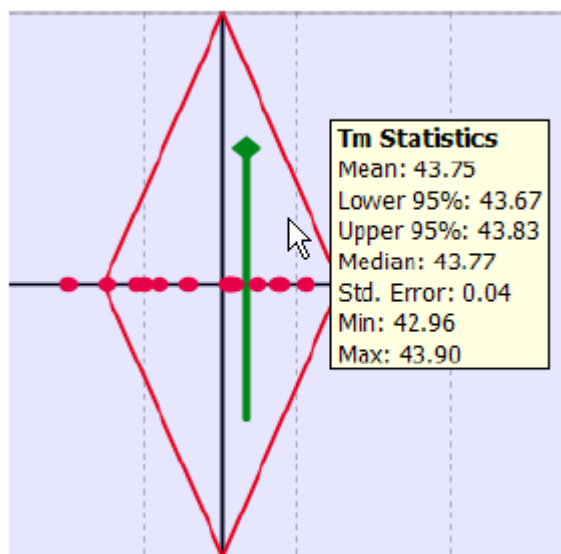
- a. Click  **Condition Hierarchy** above the top right corner of the plot.
 - b. In the dialog box, select a condition, then use the Up and Down arrows to change the hierarchy of conditions.
The condition at the top-most level of the hierarchy is displayed on the far-right side of the Replicate Results Plot, and the replicate plots are grouped according to the top-most condition.
-

Note: For the example buffer screening study, set up the hierarchy so that Buffer is at the bottom in the dialog box and displayed on the far-left side of the Replicate Results Plot.

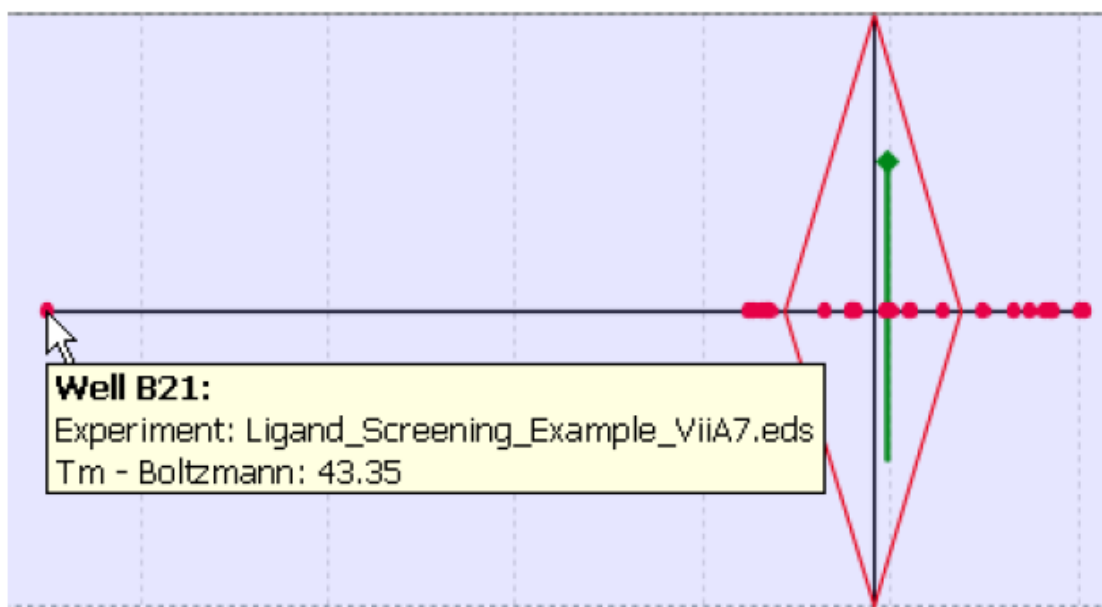
3. Scan the Replicate Results Plot to review the conditions that affect the T_m values relative to the reference replicate group.
4. Review the T_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot one or more times.
 - b. Move the plot: Click , then click-drag the plot until the replicate plot of interest is in view.




- c. Place the cursor within the diamond, then wait to view a tooltip with the T_m statistics for the replicate group.



- d. To examine outliers, place the cursor over a datapoint, then wait to view a tooltip with the well information, experiment file name, and the T_m selected for the plot.



Note: Click  to restore the default zoom.



5. In the Replicate Groups table, review the T_m statistics for each replicate group:
 - **Std. Error** (standard error of the mean for the T_m value): Is the value low? If the value is high, review the data for each replicate.
 - **Min** and **Max™** (minimum and maximum T_m values for the replicate group): Is the range of T_m values for the replicate group within 1 degree? If the range of T_m values exceeds 1 degree, review the data for each replicate.
6. Omit outliers as necessary, then click  **Analyze**.

Review the positive hits

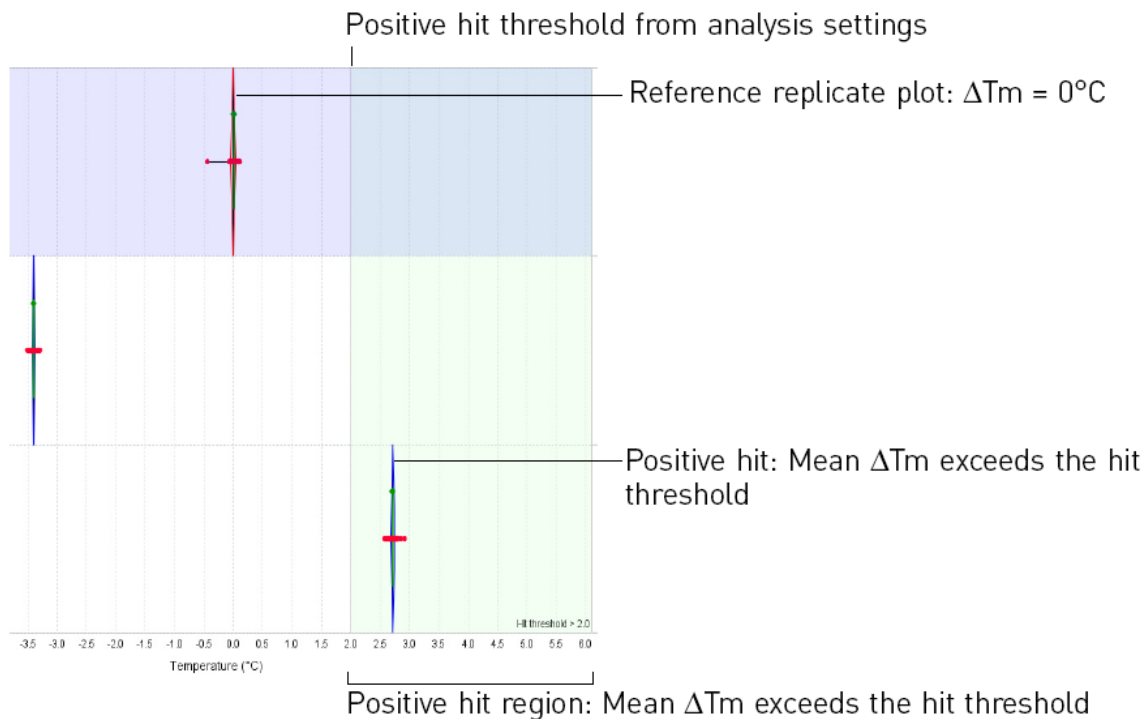
In the Analysis > Replicate Results screen, review the positive hits to identify the conditions that produce the maximum effect on thermal stability. Replicate groups with positive hits have ΔT_m values that exceed the threshold set in the analysis settings.

Note: The positive hits are determined according to the analysis settings that you specified.



Note: You must specify the reference replicate group to calculate ΔT_m values and to determine positive hits.

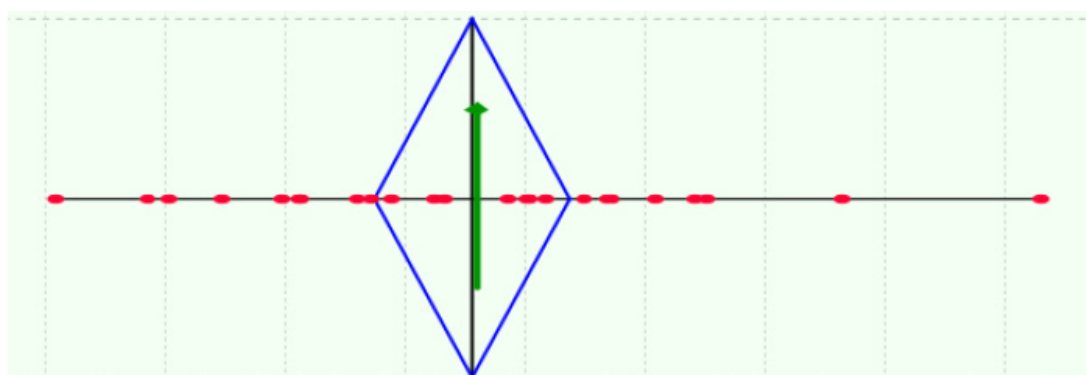
1. In the Analysis > Replicate Results screen, select  **Plot by**, then select the type of ΔT_m statistics to review:
 - **ΔT_m -Boltzmann**
 - **ΔT_m -Derivative**
2. Select the  **Show in Plot** menu, then select to show **Positive Hits**.

3. Scan the green shaded area of the Replicate Results Plot for positive hits.

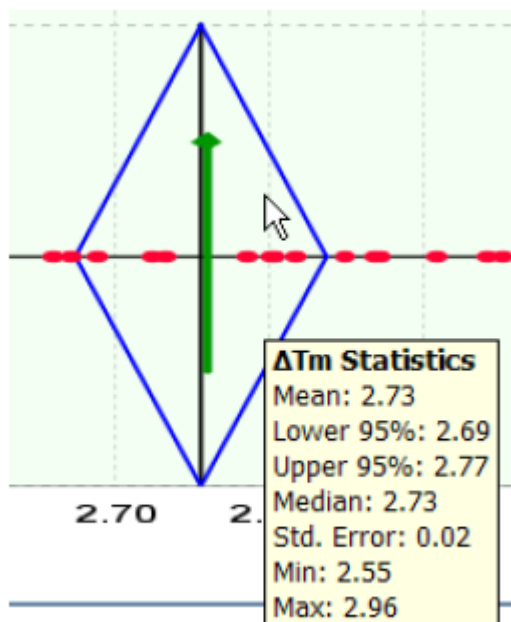


4. Review the ΔT_m statistics for each replicate group in the Replicate Results Plot:

- Zoom in: Click , then click-drag an area on the plot as many times as you need to magnify the plot.
- Move the plot: Click , then click-drag the plot until the replicate plot you want to review is in view.



- c. Place the cursor over the replicate plot, then wait to view a tooltip with the ΔT_m statistics for the replicate group.



Note: Click  to restore the default zoom.

5. In the Replicate Groups table, review the positive hits (✔) in the Hits B or Hits D column and review the ΔT_m B or ΔT_m D statistics in the table.

6

Mutation screening studies

Perform a mutation screening study to screen for mutations that increase or decrease the thermal stability of a protein.

“Create and set up an experiment file for the instrument run” on page 77



“Prepare the protein melt reactions” on page 79



“Run the protein melt reactions” on page 80



“Set up the analysis” on page 82



“Review the well results” on page 85



“Review the replicate results” on page 90

Example experiment files and plate template files are located in the examples folder: `...Program Files\Applied Biosystems\Protein Thermal Shift Software\examples`, where `...` is the installation drive.

To view the data for the buffer screening example used in this chapter, use the `Mutation_Screening_Example_ViiA7.eds` and `Mutation_Screening_Example_Setup_ViiA7.csv` files in the `ViiA7 Example Files` folder.

General guidelines

For general guidelines that apply to all Protein Thermal Shift™ studies, see “General guidelines” on page 12.

Experimental design

Before you perform a mutation screening study, we recommend that you first perform a buffer screening study to identify a buffer in which the wild type protein is thermally stable.

Replicates and controls

For buffer screening studies, we recommend that you prepare:

- At least 4 replicates of each reaction
- At least 4 replicates of no protein controls (NPCs)

Create and set up an experiment file for the instrument run

This section contains general settings for creating and setting up an experiment file. For detailed instructions, see the following sections.

Instrument	Page
QuantStudio™ 7 Pro Real-Time PCR System QuantStudio™ 6 Pro Real-Time PCR Instrument	page 15
QuantStudio™ 5 Real-Time PCR System QuantStudio™ 3 Real-Time PCR System QuantStudio™ 1 Real-Time PCR System	page 18
QuantStudio™ 7 Flex Real-Time PCR System QuantStudio™ 6 Flex Real-Time PCR System QuantStudio™ 12K Flex Real-Time PCR System	page 22
ViiA™ 7 Real-Time PCR System	page 28
StepOne™ Real-Time PCR System StepOnePlus™ Real-Time PCR System	page 31
7500 Real-Time PCR System 7500 Fast Real-Time PCR System	page 34

Experiment file settings

Setup	Setting
Experiment properties	<ul style="list-style-type: none"> • Experiment type: Melt Curve • Reagents: Other • Ramp speed: Fast or Standard
Target properties	<ul style="list-style-type: none"> • Reporter: ROX™ • Quencher: None
Plate layout	<ul style="list-style-type: none"> • Assign targets to all wells in use • Passive reference: None

(continued)

Setup	Setting
Run method	<ul style="list-style-type: none"> Reaction Volume Per Well: 20 μL Thermal profile: <ul style="list-style-type: none"> Step 1, Temp: 25°C, Time: 2 minutes Step 2, Temp: 99°C, Time: 2 minutes Ramp mode: Continuous Ramp rate: Step 1: 1.6°C/s, Step 2: 0.05°C/s for the following systems. <ul style="list-style-type: none"> QuantStudio™ 7 Pro Real-Time PCR System QuantStudio™ 6 Pro Real-Time PCR System QuantStudio™ 5 Real-Time PCR System QuantStudio™ 3 Real-Time PCR System QuantStudio™ 1 Real-Time PCR System QuantStudio™ 7 Flex Real-Time PCR System QuantStudio™ 6 Flex Real-Time PCR System QuantStudio™ 12K Flex Real-Time PCR System ViiA™ 7 Real-Time PCR System Ramp rate: 1% for the following systems. <ul style="list-style-type: none"> StepOne™ Real-Time PCR System StepOnePlus™ Real-Time PCR System 7500 Real-Time PCR System 7500 Fast Real-Time PCR System Optical Filters (ViiA™ 7 System only): <ul style="list-style-type: none"> Coupled filters: x2(520\pm10)—m2(558\pm11) Decoupled filters: x1(470\pm15)—m3(586\pm10)

Defining and assigning targets

We recommend that you define and assign targets so you can review the melt curves for the replicate groups in the Real-Time PCR System Software before importing the experiment files into the Protein Thermal Shift™ study. Define the following targets:

- One target for each buffer condition
- Define a target for the no protein control (NPC)

IMPORTANT! Make sure that the well assignments correspond exactly with the arrangement of reactions in the reaction plate. Setup errors may result in an incorrect grouping of replicates.

Example targets

In the ViiA™ 7 Real-Time PCR System mutation screening example file, targets were defined for each protein and the no protein control (NPC).

Target Name	Reporter	Quencher	Color
Wild Type	ROX	None	Red
Mutant 1	ROX	None	Blue
Mutant 2	ROX	None	Green
NPC	ROX	None	Pink

Example plate layout

In this example, columns 13–24 and rows I–P are not in use.

	1	2	3	4	5	6	7	8	9	10	11	12
A	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type	U Wild Type
B												
C	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1	U Mutant 1
D												
E	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2	U Mutant 2
F												
G												
H					U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC	U NPC

Prepare the protein melt reactions

Protein melt reaction stability

For consistency in T_m values, we recommend that you keep the protein melt reactions on ice until you are ready to load the instrument and start the run.

If the protein is thermally stable at ambient temperatures, you may consider preparing the reactions in advance and leave the reaction plate at ambient temperature, protected from light. However, the fluorescence levels will decrease over time and the T_m values will vary, depending on the protein and its thermal stability. If you want to prepare the reaction plates in advance, we recommend that you first determine the benchtop stability of your protein melt reactions.

Required materials

Required materials for protein melt reactions:

- Protein Thermal Shift™ Dye (1000X)
- Protein Thermal Shift™ Buffer
- Water
- Protein samples
- MicroAmp™ Optical Reaction Plate appropriate for your real-time PCR instrument
- MicroAmp™ Optical Adhesive Film appropriate for your reaction plate

Prepare the protein melt reactions

We recommend that you prepare four replicates of each reaction.

1. Prepare a fresh dilution of Protein Thermal Shift™ Dye (1000X) to 8X.
2. Place the appropriate reaction plate or tubes on ice, then prepare the protein melt reactions:
 - Make sure that the arrangement of reactions in the reaction plate corresponds exactly with the well assignments in the experiment file.

IMPORTANT! Setup errors may result in an incorrect grouping of the data and incorrect T_m statistics.

- Add reaction components to the plate in the order listed.

Component	Volume
Protein Thermal Shift™ Buffer	5.0 µL
Water + protein	12.5 µL
Diluted Protein Thermal Shift™ Dye (8X)	2.5 µL
Total volume for each control reaction	20.0 µL

3. Pipet each reaction up and down 10 times to mix well.
4. Seal the plate with MicroAmp™ Optical Adhesive Film, spin it at 1000 rpm for 1 minute, then place it on ice.

Run the protein melt reactions

Load and run the protein melt reactions on a supported Applied Biosystems™ Real-Time PCR System, then analyze and save the experiment file before importing the experiment file into the Protein Thermal Shift™ Software.

IMPORTANT! Keep the protein melt reactions on ice until you load the instrument.

Load and run the reactions

1. Transfer the experiment file that you created for the run to the computer that is connected to the instrument, if needed.
2. In the home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file you created for the run.
3. Load the reaction plate into the instrument.
4. In the software, start the run.

Review the melt curves

Using the Real-Time PCR System Software, open the experiment file from the completed instrument run, analyze and save the experiment file, then review the melt curves.

Note: You must analyze and save the experiment file in the Real-Time PCR System Software before you can import it into the Protein Thermal Shift™ Software.

Some common troubleshooting causes are provided here. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

1. In the Home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file from the instrument run.
2. View the melt curves:

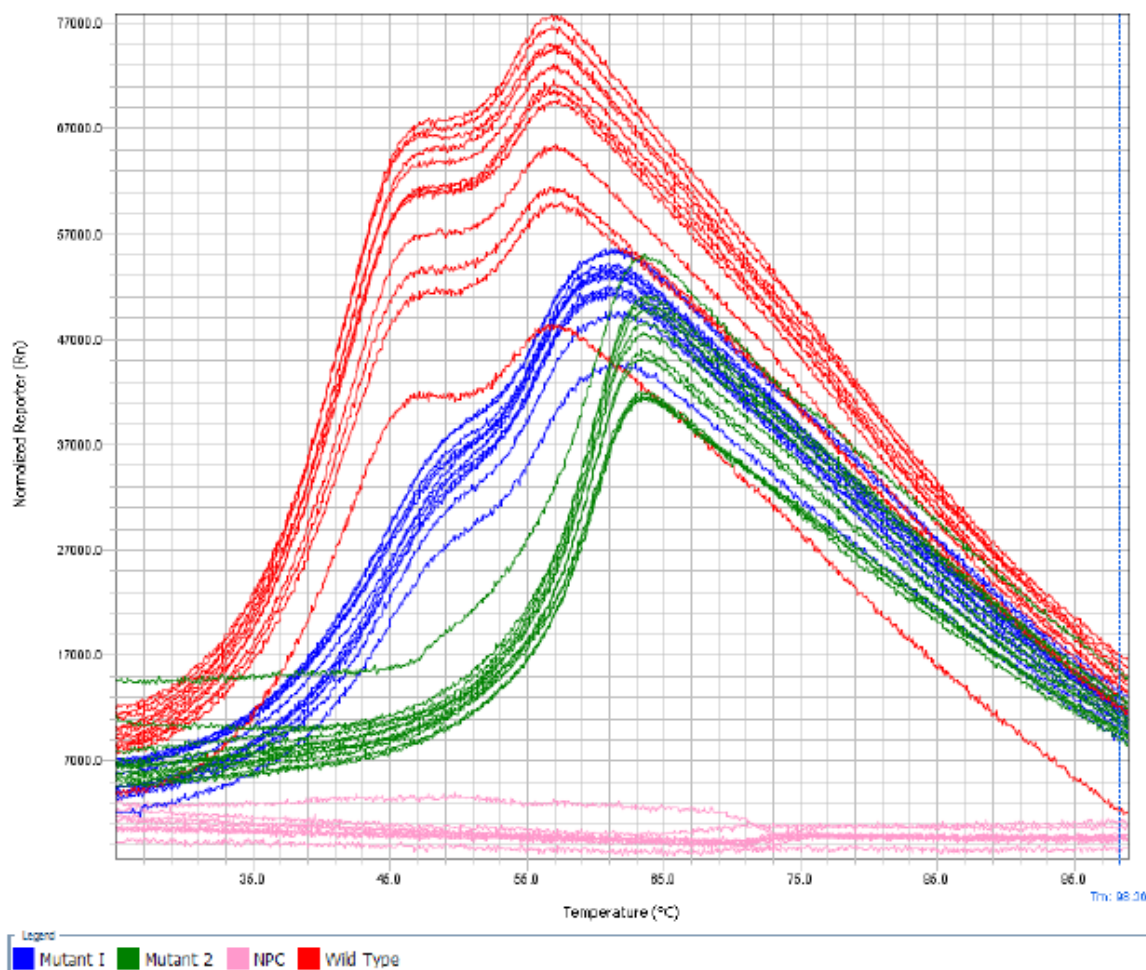
Real-time PCR System Software	View the melt curve
ViiA™ 7 Software	Click Analysis ▶ Melt Curve Plot in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.
StepOne™ Software or 7500 Software	Click Analysis ▶ Melt Curve in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.

3. Review the melt curves:
 - Do you see fluorescence signals in all of the sample wells?
No fluorescence signals in the sample wells may indicate missing dye or protein or an instrument problem.
 - Do you see flat fluorescence levels in the NPC wells?
High fluorescence levels in the NPC wells may indicate protein contamination in the wells or protein melt reactions; or it may indicate that the dye interacts with a component in the buffer.
 - Do the replicates have similar melt curves?
4. Save, then close the experiment file.

Note: The melt curves in the real-time PCR software may not exactly match the melt curves in the Protein Thermal Shift™ Software. When the experiment files are imported into the Protein Thermal Shift™ Software, the Protein Thermal Shift™ Software reduces the noise in the fluorescence data.

Example melt curves

Melt curves for the mutation screening example file from the ViiA™ 7 Real-Time PCR System:






Set up the analysis

This section provides instructions for setting up the Protein Thermal Shift™ study using Protein Thermal Shift™ Software v1.4.

Setup guidelines

- The experiment files that you import into the study must contain analyzed melt curve data from a complete melt curve run.
- Set up the analysis group so that it contains experiment files from only one instrument.

Create and set up the study



1. In the Home screen of the Protein Thermal Shift™ Software, click  **Create Study**.
2. Complete the **Setup > Properties** screen:
 - The Study Name cannot be more than 100 characters and cannot contain these characters: / \ * " ? < > | . ,
 - The instrument selection must match the instrument type that you used to run the protein melt reactions and generate the experiment files.
3. In the **Setup > Conditions** screen, define the conditions and condition values, then list the analysis groups for your study.
 - a. In the **Condition Names** pane, click  **Add** to add up to 20 custom parameters for the study, such as pH.
 - b. Select a condition of interest, then in the pane for the values for the condition, add values for the condition.
 - c. In the **Analysis Groups** pane, create groups, then assign wells to each group to perform a more focused analysis with the software, if needed.
4. In the **Setup > Experiment Files** screen, click  **Import**, then select the experiment file (EDS) for the instrument type that you selected for the study. Repeat for each experiment file to import into the study.

Note: The experiment file name and data are imported. Plate setup information is not imported.

5. For each experiment file, assign the conditions, task, and analysis group to the wells that contained protein melt reactions.
 - a. In the **Setup > Experiment Files** screen, click-drag wells in the plate to select multiple wells. If the wells are not continuous, ctrl+click the wells of interest.
 - b. Right-click one of the selected wells, then click **Assign Well Content**.
 - c. In the dialog box that appears, select the value for each condition that applies to the selected wells, then click **Done**.
In the **Task** dropdown list, you can select **Sample**, **Reference** or **NPC** to assign the well as an unknown sample, a reference sample, or a no protein control.
 - d. Repeat step 5 and substep 5a for each condition.
You can click **Auto-Fill Settings** to assign the wells faster.

6. Ensure that the condition assignments correspond exactly with the contents of the reaction plate.

IMPORTANT! Setup errors can result in an incorrect grouping of the data and incorrect T_m statistics.

7. For each analysis group, assign the reference replicate group:
 - a. Select the wells for the replicate group to use as the reference.
 - b. Click  **Assign**, then select **Reference** as the Task.
8. Click  **Save** in the toolbar to save and analyze the study.


Assign wells automatically

You can use auto fill settings to assign wells faster.


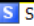

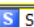
Alternatively, you can click **Load Plate Template** to import a Microsoft™ Excel™ spreadsheet with the well information.

Create and set up the study.

1. In the **Setup > Experiment Files** screen, assign one or more conditions to one or more rows or columns of wells.
2. Click **Auto-Fill Settings ▶ Dock**.
3. In the **Auto-Fill Settings** pane, make the following selections for each condition.
 - a. Click the ☒ (**checkmark**) to deselect a condition so that values are not automatically assigned for the condition, if needed.
 - b. For each selected condition, select **Copy** to copy the value for the condition, or select **Series** to extend the sequence for the series of values for the condition.

For each series, the values are added in the order in which the values were created in the study setup.
4. For each column or row that you want to fill, click-drag to select wells from which the condition values are copied or extended.
5. Place the pointer over the last selected well of the column or row, then drag the pointer over the wells that you want to fill.
6. Click  **Save** to save the study.

The remaining wells are filled according to the selections.


7	8	9	10
 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1		
 Sample Protein A 100 mM AG 1	 Sample Protein A 150 mM AG 1		

Well 18
Task: Sample
Analysis Group: AG 1
Protein: Protein A
Salt: 150 mM

Review the analysis settings

Review and edit the analysis settings to optimize the analysis for your study.

For the examples shown in this user guide, no changes were made to the analysis settings. If you are reviewing the data in the example studies provided with the software, try revising the analysis settings to see how the settings affect the positive hits and the flags.

1. Click  **Analysis Settings** in the toolbar.
2. On the Positive Hit tab, specify the ΔT_m -Boltzmann and ΔT_m -Derivative values to indicate a positive hit:
 - Select **>** to identify buffer conditions that increase protein thermal stability, or select **<** to identify buffer conditions that decrease protein thermal stability.
 - Enter the number of degrees ($^{\circ}\text{C}$) of T_m shift relative to the reference to indicate a positive hit.
3. On the Flags tab, specify settings for applying flags:
 - a. Select the flags to use in the analysis.
 - b. For the High Background, High NPC, Low Signal, and Poor Fit flags, specify the condition and threshold for applying the flag.
4. Apply the analysis settings and analyze:
 - Click **Apply** to apply the analysis settings and reanalyze while keeping the Analysis Settings dialog box open.
or
 - Click **OK** to apply the analysis settings, reanalyze, and close the Analysis Settings.

Review the well results

Using the Protein Thermal Shift™ Software, review the melt curves and well table and optimize the analysis in the Analysis > Well Results screen.



The Well Results screen displays fluorescence and derivative melt curve plots, calculated T_m values, individual well results, and flags. As necessary, edit the analysis settings, edit the baseline, edit the region of analysis, edit the analysis mode, and omit outliers.

This section provides guidance on how to review and interpret the well results. Some common troubleshooting causes are provided. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

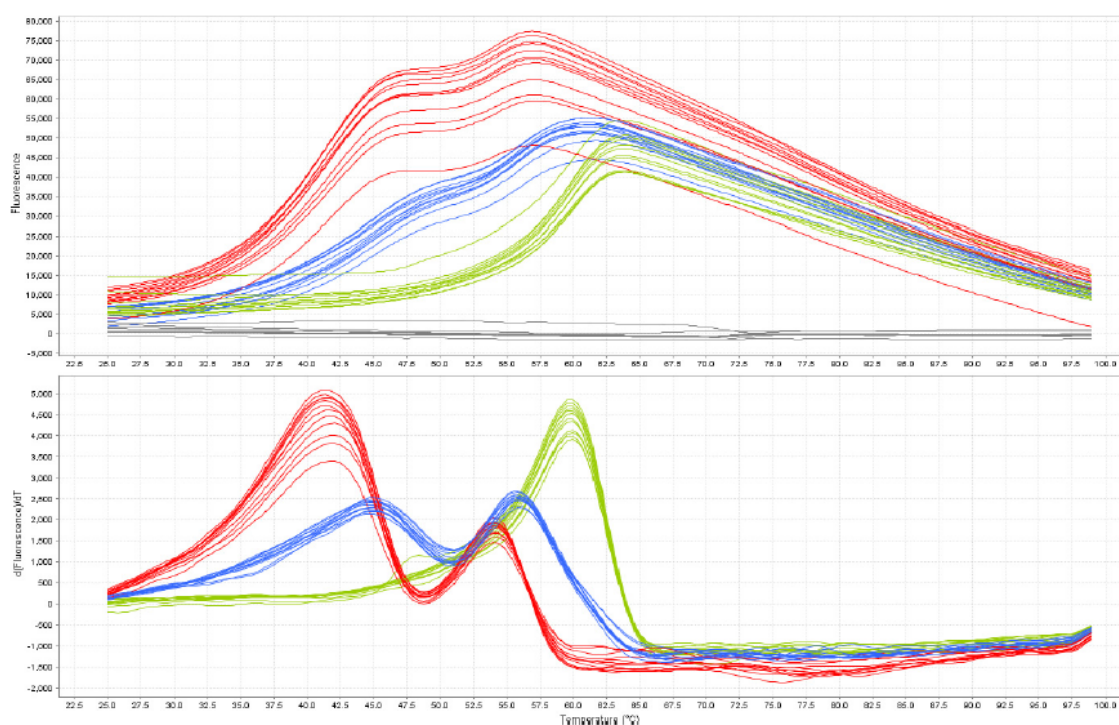
Review the melt curves



Review the melt curve plots to visualize the fluorescence and derivative fluorescence data. If necessary, change the analysis mode.

Note: For NPC wells, the derivative melt curves are not displayed.

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Unselected Wells
 - Legend
2. Select  **Color By**, then select **Protein** to color the melt curves according to the protein condition value assigned for the well.

Note: To set the color for each protein variant to make the plots easier to distinguish, go to the Setup > Conditions screen.




3. For each replicate group, select all the wells in the replicate group, then review the fluorescence levels in the melt curves:
 - For the NPC wells, do you observe a rise in fluorescence? If so, the wells or protein melt reactions may be contaminated with protein or the dye may interact with a buffer component.
 - For sample or reference wells, do you observe flat melt curves? If so, condition assignments may be incorrect or a component is missing from the protein melt reactions.
 - Within each replicate group, are the fluorescence melt curves similar to each other? Within each replicate group, are the derivative melt curves similar to each other? If the melt curves for the replicates are dissimilar, pipetting errors may have occurred during reaction setup or condition assignments may be incorrect.
4. If the derivative melt curves for the replicate group show multiple melt phases, set the analysis mode to Auto: Multiple T_m, then review the derivative melt curves:
 - a. In the Well Table or in the melt curve plot, select the wells with multiple melt phases.
 - b. Click  **Auto Analysis Options**, then select **Auto: Multiple T_m**.
 - c. Click  **Analyze** to reanalyze using the Auto: Multiple T_m analysis mode.
 - d. Review the number of melt phases in the derivative melt curves:
 - Do all replicate groups have the same number of melt phases?
 - For each replicate group, are there outliers with a different number of melt phases than the other samples in the replicate group? You may consider omitting outliers from analysis.

Review the regions of analysis (ROA)

Review the ROAs detected by the software. If necessary, edit the ROAs.

Note: If no melt phases are detected, no ROAs are defined. Negative controls should have no melt phases and no ROAs.

1. In the Analysis > Well Results screen, confirm that each ROA meets the following criteria:
 - For melt curves with one melt phase, the curve within the ROA resembles a sigmoidal profile.
 - At the start temperature, the signal is relatively flat.
 - At the end temperature, the signal has already reached its maximum.
2. For each replicate group, edit the ROAs so that all of the wells in the replicate group have the same ROAs:
 - a. Select the replicates, Click  (Define ROA) in the toolbar above the melt curve plots, then click-drag an area in one of the plots to define a melt phase and replace the ROA. Repeat for each melt phase you identify.

- b. If necessary, adjust the start and end temperatures for each ROA:
 - **To move the ROA:** Starting from within the ROA, click-drag the ROA.
 - **To adjust the start and end temperatures individually:** Click-drag the ROA start or the end line.



IMPORTANT! Make sure that the fluorescence at the start temperature is lower than the fluorescence at the end temperature.

3. Click ► **Analyze** to reanalyze using the edited ROAs.

Note: After you edit ROAs, the analysis mode changes from Auto to Manual.

Review the flags and T_m values

Review the flags and T_m values, then consider editing the analysis settings and/or omitting wells before you review the replicate results.

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann T_m
 - Derivative T_m
 - Unselected Wells
 - Legend
2. Review flagged wells in the Well Table:
 - a. Click the  (Flag Indicator) column header to sort the wells according to the number of flags applied to the well.
 - b. Scroll the table to the right to view the flags that are applied to the wells.




- c. For flagged wells, select the well in the Well Table, then review the melt curves for the well, compared to the other wells in the replicate group.
- d. Omit wells from analysis as necessary.

3. For each replicate group, select all the wells in the replicate group, then review the Tm B (Boltzmann Tm) and the Tm D (derivative Tm) in the Well Table and in the melt curves:

Note: In the melt curves, the Boltzmann Tm is a green dashed vertical line, and the derivative Tm is black dotted vertical line.

- Are the Tm B or Tm D values significantly different from the Tm values for other wells in the replicate group? Do any replicates have melt curves that are inconsistent with the other melt curves for the replicate group? If so, you may consider omitting wells from analysis.
 - Are the melt curves within the replicate group similar?
4. If you omitted any wells from the analysis, click ► **Analyze**.

Review the Boltzmann fit

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann fit
 - Unselected Wells
 - Legend
2. Scroll through each well in the Well Table, then compare the fluorescence melt curve to the Boltzmann fit curve (dark green thick curve) and review the value in the B Fit (Boltzmann fit) column of the Well Table:
 - How well does the fluorescence melt curve correspond to the Boltzmann fit curve?
 - Is the B Fit value close to 1?
 - Is the B Fit value similar among wells in the replicate group?

Note: When you define the ROA manually, you may observe a gap between the ROA start or end temperature and the start or end of the Boltzmann fit curve. The gap occurs if the defined ROA start or end temperature does not correspond exactly with a fluorescence datapoint.

Example well results

For the ViiA™ 7 Real-Time PCR System example, observe the following:

- Fluorescence levels are flat in the NPC wells.
- For the sample and reference replicate groups, the fluorescence levels as displayed in the fluorescence melt curve are not significantly different, so you do not need to edit the baseline.
- The sample and reference wells contain one peak in the derivative melt curve.
- Within each replicate group:
 - The fluorescence melt curves and the Boltzmann Tm values are similar.
 - The derivative melt curves and the derivative Tm values are similar and there are no outliers.
- For each well, the region of analysis meets the recommended criteria.


Review the replicate results

Using the Protein Thermal Shift™ Software, review the T_m statistics for the replicate groups and look for positive hits in the Analysis > Replicate Results screen.


This section provides guidance on how to review and interpret the replicate results. Some common troubleshooting causes are provided. For more detailed information on troubleshooting, see Chapter 8, “Troubleshooting”.

Review the T_m statistics



In the Analysis > Replicate Results screen, review the T_m statistics to evaluate the variability among replicates. The T_m statistics for each ROA are plotted along the x-axis for each replicate group.

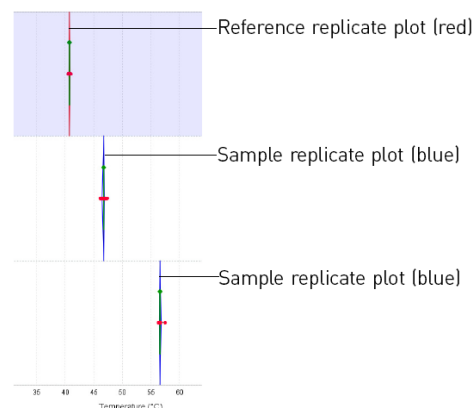
1. In the Analysis > Replicate Results screen, select  **Plot by**, then select the type of T_m statistics to review:
 - **T_m-Boltzmann**
 - **T_m-Derivative**
2. Specify the condition hierarchy to group the replicate plots and change the order of conditions in the plot:

Note: Changing the condition hierarchy does not affect the results; it only changes how the replicate plots are grouped and the order in which the conditions are displayed in the Replicate Results Plot.

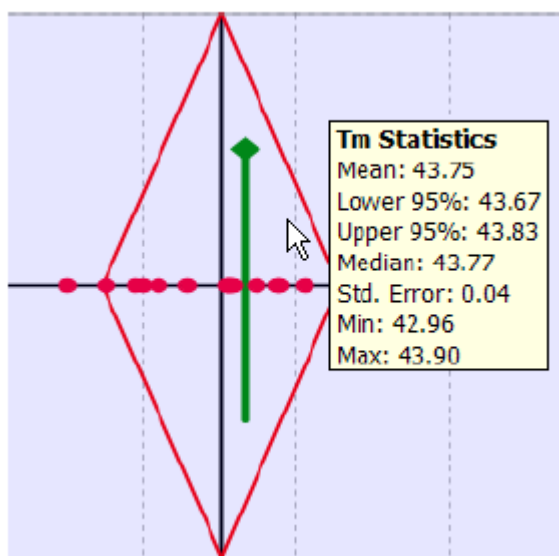
- a. Click  **Condition Hierarchy** above the top right corner of the plot.
- b. In the dialog box, select a condition, then use the Up and Down arrows to change the hierarchy of conditions.
The condition at the top-most level of the hierarchy is displayed on the far-right side of the Replicate Results Plot, and the replicate plots are grouped according to the top-most condition.

Note: For the example mutation screening study, set up the hierarchy so that Protein is at the bottom in the dialog box and displayed on the far-left side of the Replicate Results Plot.

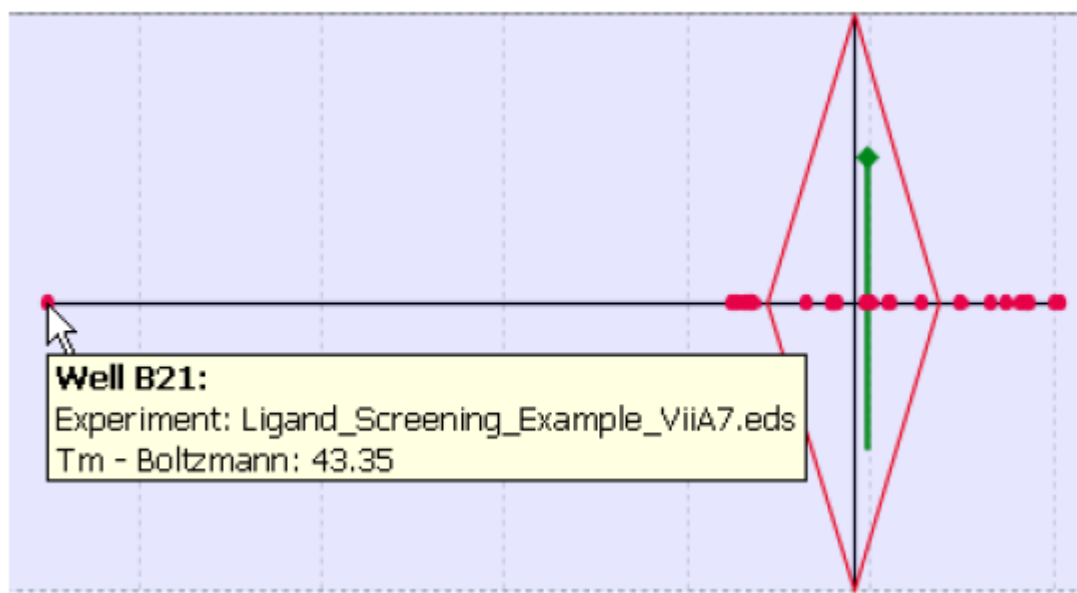
3. Scan the Replicate Results Plot to review the conditions that affect the T_m values relative to the reference replicate group.
4. Review the T_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot one or more times.
 - b. Move the plot: Click , then click-drag the plot until the replicate plot of interest is in view.



- c. Place the cursor within the diamond, then wait to view a tooltip with the T_m statistics for the replicate group.



- d. To examine outliers, place the cursor over a datapoint, then wait to view a tooltip with the well information, experiment file name, and the T_m selected for the plot.



Note: Click to restore the default zoom. [🔍](#)



5. In the Replicate Groups table, review the T_m statistics for each replicate group:
 - **Std. Error** (standard error of the mean for the T_m value): Is the value low? If the value is high, review the data for each replicate.
 - **Min** and **Max**[™] (minimum and maximum T_m values for the replicate group): Is the range of T_m values for the replicate group within 1 degree? If the range of T_m values exceeds 1 degree, review the data for each replicate.
6. Omit outliers as necessary, then Click ► **Analyze**.

Review the positive hits

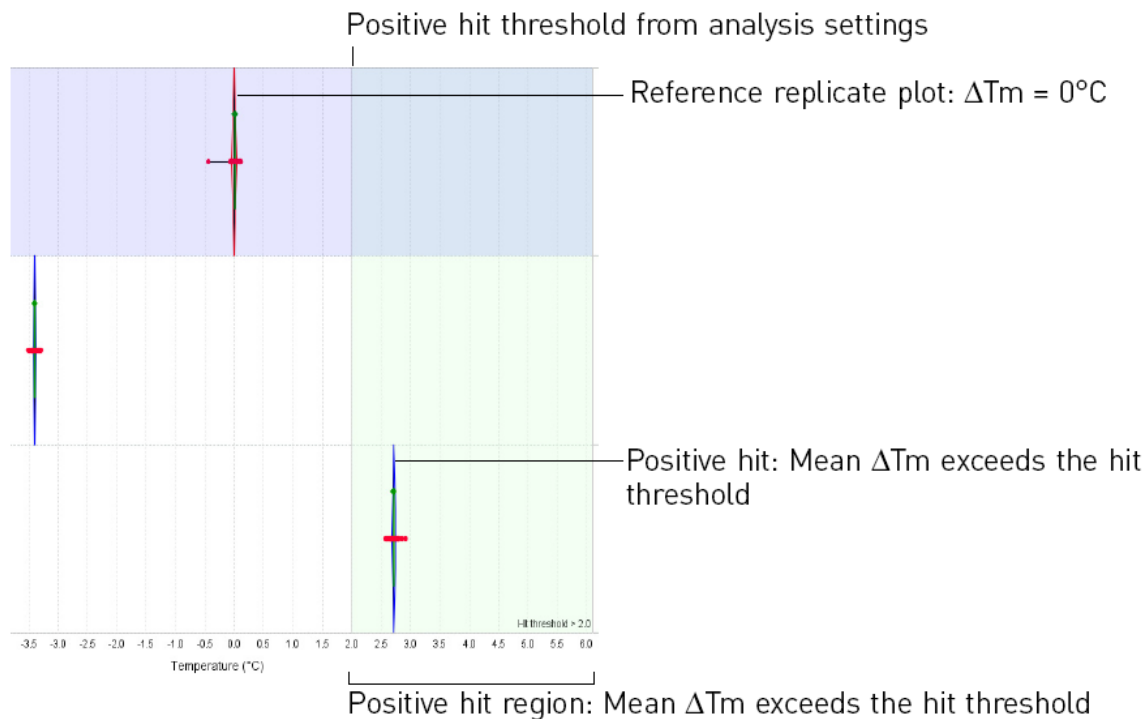
In the Analysis > Replicate Results screen, review the positive hits to identify the conditions that produce the maximum effect on thermal stability. Replicate groups with positive hits have ΔT_m values that exceed the threshold set in the analysis settings.



Note: The positive hits are determined according to the analysis settings that you specified.

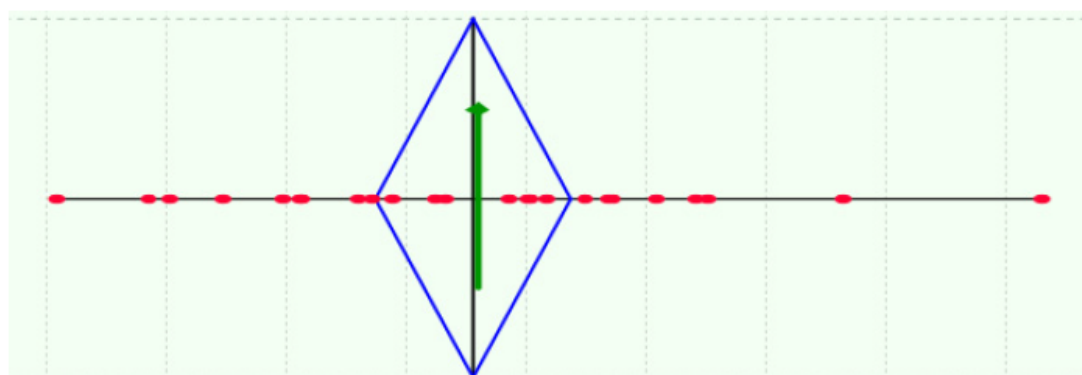
Note: You must specify the reference replicate group to calculate ΔT_m values and to determine positive hits.

1. In the Analysis > Replicate Results screen, select  **Plot by**, then select the type of ΔT_m statistics to review:
 - **ΔT_m -Boltzmann**
 - **ΔT_m -Derivative**
2. Select the  **Show in Plot** menu, then select to show **Positive Hits**.

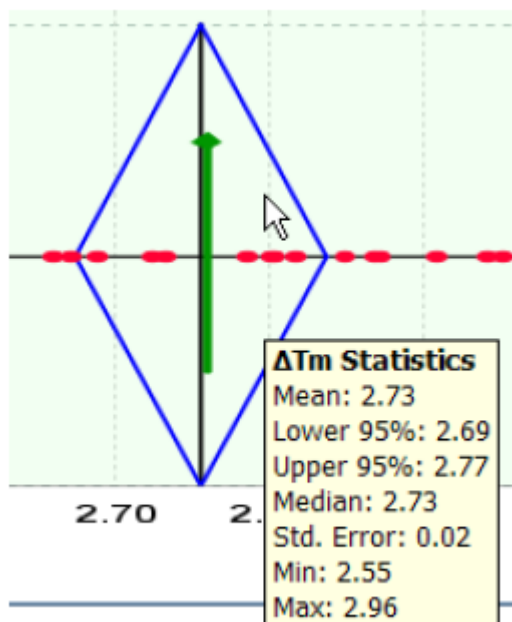
3. Scan the green shaded area of the Replicate Results Plot for positive hits.



4. Review the ΔT_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot as many times as you need to magnify the plot.
 - b. Move the plot: Click , then click-drag the plot until the replicate plot you want to review is in view.



- c. Place the cursor over the replicate plot, then wait to view a tooltip with the ΔT_m statistics for the replicate group.



Note: Click  to restore the default zoom.

5. In the Replicate Groups table, review the positive hits (🟢) in the Hits B or Hits D column and review the ΔT_m B or ΔT_m D statistics in the table.

7

Ligand screening studies

Perform a ligand screening study to identify ligand candidates that stabilize a protein and lead to successful protein crystallization.

“Create and set up an experiment file for the instrument run” on page 96



“Prepare the protein melt reactions” on page 98



“Run the protein melt reactions” on page 99



“Set up the analysis” on page 101



“Review the well results” on page 104



“Review the replicate results” on page 109

Example experiment files and plate template files are located in the examples folder: `...Program Files\Applied Biosystems\Protein Thermal Shift Software\examples`, where `...` is the installation drive.

To view the data for the ligand screening example used in this chapter, use the `Ligand_Screening_Example_ViiA7.eds` and `Ligand_Screening_Example_Setup_ViiA7.csv` files in the `ViiA7 Example Files` folder.

General guidelines

For general guidelines that apply to all Protein Thermal Shift™ studies, see “General guidelines” on page 12.

Experimental design

Before you perform a ligand screening study, we recommend that you first perform a buffer screening study to identify a buffer in which the protein is thermally stable.

After you identify ligand candidates, you can perform a ligand titration study to determine the optimal ligand:protein ratio.

Replicates and controls

For ligand screening studies, we recommend that you prepare:

- At least 4 replicates of each reaction
- At least 4 replicates of no protein controls (NPCs)
- At least 4 replicates of ligand only controls (LOCs)

Create and set up an experiment file for the instrument run

This section contains general settings for creating and setting up an experiment file. For detailed instructions, see the following sections.

Instrument	Page
QuantStudio™ 7 Pro Real-Time PCR System QuantStudio™ 6 Pro Real-Time PCR Instrument	page 15
QuantStudio™ 5 Real-Time PCR System QuantStudio™ 3 Real-Time PCR System QuantStudio™ 1 Real-Time PCR System	page 18
QuantStudio™ 7 Flex Real-Time PCR System QuantStudio™ 6 Flex Real-Time PCR System QuantStudio™ 12K Flex Real-Time PCR System	page 22
ViiA™ 7 Real-Time PCR System	page 28
StepOne™ Real-Time PCR System StepOnePlus™ Real-Time PCR System	page 31
7500 Real-Time PCR System 7500 Fast Real-Time PCR System	page 34

Experiment file settings

Setup	Setting
Experiment properties	<ul style="list-style-type: none"> • Experiment type: Melt Curve • Reagents: Other • Ramp speed: Fast or Standard
Target properties	<ul style="list-style-type: none"> • Reporter: ROX™ • Quencher: None
Plate layout	<ul style="list-style-type: none"> • Assign targets to all wells in use • Passive reference: None

(continued)

Setup	Setting
Run method	<ul style="list-style-type: none"> Reaction Volume Per Well: 20 μL Thermal profile: <ul style="list-style-type: none"> Step 1, Temp: 25°C, Time: 2 minutes Step 2, Temp: 99°C, Time: 2 minutes Ramp mode: Continuous Ramp rate: Step 1: 1.6°C/s, Step 2: 0.05°C/s for the following systems. <ul style="list-style-type: none"> QuantStudio™ 7 Pro Real-Time PCR System QuantStudio™ 6 Pro Real-Time PCR System QuantStudio™ 5 Real-Time PCR System QuantStudio™ 3 Real-Time PCR System QuantStudio™ 1 Real-Time PCR System QuantStudio™ 7 Flex Real-Time PCR System QuantStudio™ 6 Flex Real-Time PCR System QuantStudio™ 12K Flex Real-Time PCR System ViiA™ 7 Real-Time PCR System Ramp rate: 1% for the following systems. <ul style="list-style-type: none"> StepOne™ Real-Time PCR System StepOnePlus™ Real-Time PCR System 7500 Real-Time PCR System 7500 Fast Real-Time PCR System Optical Filters (ViiA™ 7 System only): <ul style="list-style-type: none"> Coupled filters: x2(520\pm10)—m2(558\pm11) Decoupled filters: x1(470\pm15)—m3(586\pm10)

Defining and assigning targets

We recommend that you define and assign targets so you can review the melt curves for the replicate groups in the Real-Time PCR System Software before importing the experiment files into the Protein Thermal Shift™ study. Define the following targets:

- One target for each ligand or for each ligand titration
- Define a target for the no protein control (NPC)
- Define a target for the ligand only control (LOC)

IMPORTANT! Make sure that the well assignments correspond exactly with the arrangement of reactions in the reaction plate. Setup errors may result in an incorrect grouping of replicates.

Example targets

In the ViiA™ 7 Real-Time PCR System ligand screening example file, targets were defined for each concentration of ligand, the ligand only control (LOC), and no protein control (NPC).

Target Name	Reporter	Quencher	Color
ProteinA +0 mM L	ROX	None	Green
ProteinA +0.1 mM L	ROX	None	Red
ProteinA +1 mM L	ROX	None	Blue
LOC	ROX	None	Pink
NPC	ROX	None	Grey

Example plate layout

In this example, rows F–P are not in use.

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L	ProteinA +0 mM L
B	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L	ProteinA +0.1 mM L
C	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L	ProteinA +1 mM L
D	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC	LOC
E	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC	NPC

Prepare the protein melt reactions

Protein melt reaction stability

For consistency in T_m values, we recommend that you keep the protein melt reactions on ice until you are ready to load the instrument and start the run.

If the protein is thermally stable at ambient temperatures, you may consider preparing the reactions in advance and leave the reaction plate at ambient temperature, protected from light. However, the fluorescence levels will decrease over time and the T_m values will vary, depending on the protein and its thermal stability. If you want to prepare the reaction plates in advance, we recommend that you first determine the benchtop stability of your protein melt reactions.

Required materials

Required materials for protein melt reactions:

- Protein Thermal Shift™ Dye (1000X)
- Protein Thermal Shift™ Buffer
- Water
- Protein
- Ligand titrations

- MicroAmp™ Optical Reaction Plate appropriate for your real-time PCR instrument
- MicroAmp™ Optical Adhesive Film appropriate for your reaction plate

Prepare the protein melt reactions

We recommend that you prepare four replicates of each reaction.

1. Prepare a fresh dilution of DyeProtein Thermal Shift™ Dye (1000X) to 8X.
2. Place the appropriate reaction plate or tubes on ice, then prepare the protein melt reactions:
 - Make sure that the arrangement of reactions in the reaction plate corresponds exactly with the well assignments in the experiment file.

IMPORTANT! Setup errors may result in an incorrect grouping of the data and incorrect T_m statistics.

- Add reaction components to the plate in the order listed.

Component	Volume
Protein Thermal Shift™ Buffer	5.0 µL
Water + protein + ligand	12.5 µL
Diluted Protein Thermal Shift™ Dye (8X)	2.5 µL
Total volume for each control reaction	20.0 µL

3. Pipet each reaction up and down 10 times to mix well.
4. Seal the plate with MicroAmp™ Optical Adhesive Film, spin it at 1000 rpm for 1 minute, then place it on ice.

Run the protein melt reactions

Load and run the protein melt reactions on a supported Applied Biosystems™ Real-Time PCR System, then analyze and save the experiment file before importing the experiment file into the Protein Thermal Shift™ Software.

IMPORTANT! Keep the protein melt reactions on ice until you load the instrument.

Load and run the reactions

1. Transfer the experiment file that you created for the run to the computer that is connected to the instrument, if needed.
2. In the home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file you created for the run.
3. Load the reaction plate into the instrument.
4. In the software, start the run.

Review the melt curves

Using the Real-Time PCR System Software, open the experiment file from the completed instrument run, analyze and save the experiment file, then review the melt curves.

Note: You must analyze and save the experiment file in the Real-Time PCR System Software before you can import it into the Protein Thermal Shift™ Software.

Some common troubleshooting causes are provided here. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

1. In the Home screen of the Real-Time PCR System Software, click **Open**, then select the experiment file from the instrument run.
2. View the melt curves:

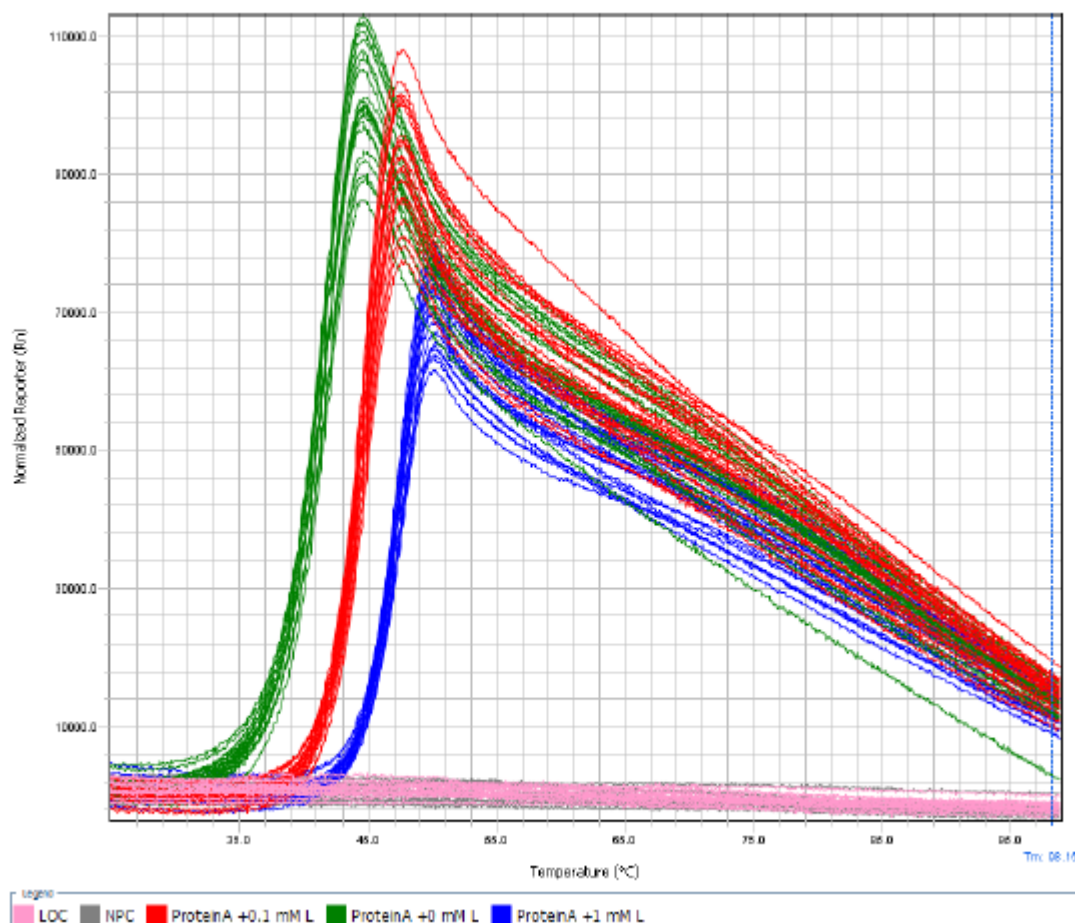
Real-time PCR System Software	View the melt curve
ViiA™ 7 Software	Click Analysis ▶ Melt Curve Plot in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.
StepOne™ Software or 7500 Software	Click Analysis ▶ Melt Curve in the navigation pane, select Normalized Reporter from the Plot dropdown list, then select Target from the Color dropdown list.

3. Review the melt curves:
 - Do you see fluorescence signals in all of the sample wells?
No fluorescence signals in the sample wells may indicate missing dye or protein or an instrument problem.
 - Do you see flat fluorescence levels in the NPC wells?
High fluorescence levels in the NPC wells may indicate protein contamination in the wells or protein melt reactions; or it may indicate that the dye interacts with a component in the buffer.
 - Do the replicates have similar melt curves?
4. Save, then close the experiment file.

Note: The melt curves in the real-time PCR software may not exactly match the melt curves in the Protein Thermal Shift™ Software. When the experiment files are imported into the Protein Thermal Shift™ Software, the Protein Thermal Shift™ Software reduces the noise in the fluorescence data.

Example melt curves

Melt curves for the ligand screening example file from the ViiA™ 7 Real-Time PCR System:






Set up the analysis

This section provides instructions for setting up the Protein Thermal Shift™ study using Protein Thermal Shift™ Software v1.4.

Setup guidelines

- The experiment files that you import into the study must contain analyzed melt curve data from a complete melt curve run.
- Set up the analysis group so that it contains experiment files from only one instrument.

Create and set up the study

1. In the Home screen of the Protein Thermal Shift™ Software, click  **Create Study**.
2. Complete the **Setup > Properties** screen:
 - The Study Name cannot be more than 100 characters and cannot contain these characters: / \ * “ ? < > | . ,
 - The instrument selection must match the instrument type that you used to run the protein melt reactions and generate the experiment files.
3. In the **Setup > Conditions** screen, define the conditions and condition values, then list the analysis groups for your study.
 - a. In the **Condition Names** pane, click  **Add** to add up to 20 custom parameters for the study, such as pH.
 - b. Select a condition of interest, then in the pane for the values for the condition, add values for the condition.
 - c. In the **Analysis Groups** pane, create groups, then assign wells to each group to perform a more focused analysis with the software, if needed.
4. In the **Setup > Experiment Files** screen, click  **Import**, then select the experiment file (EDS) for the instrument type that you selected for the study. Repeat for each experiment file to import into the study.

Note: The experiment file name and data are imported. Plate setup information is not imported.

5. For each experiment file, assign the conditions, task, and analysis group to the wells that contained protein melt reactions.
 - a. In the **Setup > Experiment Files** screen, click-drag wells in the plate to select multiple wells. If the wells are not continuous, ctrl+click the wells of interest.
 - b. Right-click one of the selected wells, then click **Assign Well Content**.
 - c. In the dialog box that appears, select the value for each condition that applies to the selected wells, then click **Done**.
In the **Task** dropdown list, you can select **Sample**, **Reference** or **NPC** to assign the well as an unknown sample, a reference sample, or a no protein control.
 - d. Repeat step 5 and substep 5a for each condition.
You can click **Auto-Fill Settings** to assign the wells faster.
6. Ensure that the condition assignments correspond exactly with the contents of the reaction plate.


IMPORTANT! Setup errors can result in an incorrect grouping of the data and incorrect T_m statistics.

7. For each analysis group, assign the reference replicate group:
 - a. Select the wells for the replicate group to use as the reference.

Review the analysis settings

Review and edit the analysis settings to optimize the analysis for your study.

For the examples shown in this user guide, no changes were made to the analysis settings. If you are reviewing the data in the example studies provided with the software, try revising the analysis settings to see how the settings affect the positive hits and the flags.

1. Click  **Analysis Settings** in the toolbar.
2. On the Positive Hit tab, specify the ΔT_m -Boltzmann and ΔT_m -Derivative values to indicate a positive hit:
 - Select **>** to identify buffer conditions that increase protein thermal stability, or select **<** to identify buffer conditions that decrease protein thermal stability.
 - Enter the number of degrees ($^{\circ}\text{C}$) of T_m shift relative to the reference to indicate a positive hit.
3. On the Flags tab, specify settings for applying flags:
 - a. Select the flags to use in the analysis.
 - b. For the High Background, High NPC, Low Signal, and Poor Fit flags, specify the condition and threshold for applying the flag.
4. Apply the analysis settings and analyze:
 - Click **Apply** to apply the analysis settings and reanalyze while keeping the Analysis Settings dialog box open.
or
 - Click **OK** to apply the analysis settings, reanalyze, and close the Analysis Settings.

Review the well results

Using the Protein Thermal Shift™ Software, review the melt curves and well table and optimize the analysis in the Analysis > Well Results screen.



The Well Results screen displays fluorescence and derivative melt curve plots, calculated T_m values, individual well results, and flags. As necessary, edit the analysis settings, edit the baseline, edit the region of analysis, edit the analysis mode, and omit outliers.

This section provides guidance on how to review and interpret the well results. Some common troubleshooting causes are provided. For more detailed troubleshooting information, see Chapter 8, “Troubleshooting”.

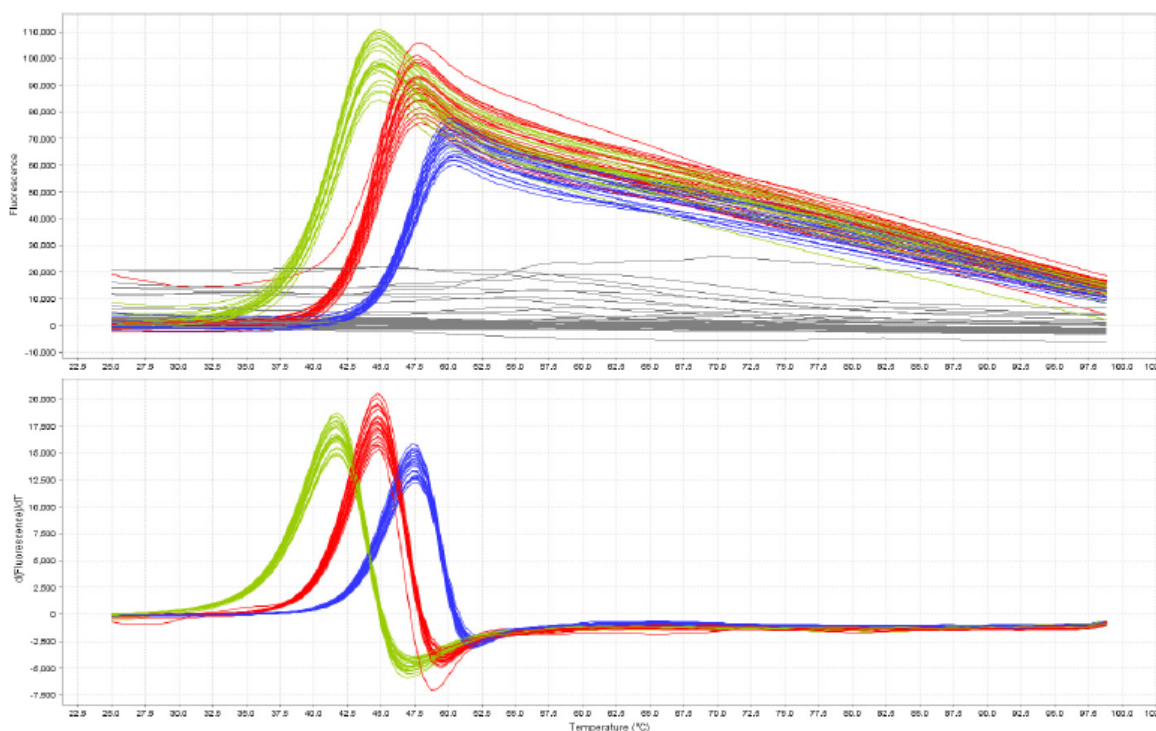
Review the melt curves



Review the melt curve plots to visualize the fluorescence and derivative fluorescence data. If necessary, change the analysis mode.

Note: For NPC wells, the derivative melt curves are not displayed.

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Unselected Wells
 - Legend
2. Select  **Color By**, then select **Ligand** to color the melt curves according to the ligand condition value assigned for the well.

Note: To set the color for each ligand to make the plots easier to distinguish, go to the Setup > Conditions screen.




3. For each replicate group, select all the wells in the replicate group, then review the fluorescence levels in the melt curves:
 - For the NPC wells, do you observe a rise in fluorescence? If so, the wells or protein melt reactions may be contaminated with protein or the dye may interact with a buffer component.
 - For the LOC wells, do you observe a rise in fluorescence? A rise in fluorescence in LOC wells but not in NPC wells may indicate protein contamination in the ligand or ligand-dye interactions.
 - For sample or reference wells, do you observe flat melt curves? If so, condition assignments may be incorrect or a component is missing from the protein melt reactions.
 - Within each replicate group, are the fluorescence melt curves similar to each other? Within each replicate group, are the derivative melt curves similar to each other? If the melt curves for the replicates are dissimilar, pipetting errors may have occurred during reaction setup or condition assignments may be incorrect.
4. If the derivative melt curves for the replicate group show multiple melt phases, set the analysis mode to Auto: Multiple T_m, then review the derivative melt curves:
 - a. In the Well Table or in the melt curve plot, select the wells with multiple melt phases.
 - b. Click  **Auto Analysis Options**, then select **Auto: Multiple T_m**.
 - c. Click  **Analyze** to reanalyze using the Auto: Multiple T_m analysis mode.
 - d. Review the number of melt phases in the derivative melt curves:
 - Do all replicate groups have the same number of melt phases?
 - For each replicate group, are there outliers with a different number of melt phases than the other samples in the replicate group? You may consider omitting outliers from analysis.

Review the regions of analysis (ROA)

Review the ROAs detected by the software. If necessary, edit the ROAs.

Note: If no melt phases are detected, no ROAs are defined. Negative controls should have no melt phases and no ROAs.

1. In the Analysis > Well Results screen, confirm that each ROA meets the following criteria:
 - For melt curves with one melt phase, the curve within the ROA resembles a sigmoidal profile.
 - At the start temperature, the signal is relatively flat.
 - At the end temperature, the signal has already reached its maximum.
2. For each replicate group, edit the ROAs so that all of the wells in the replicate group have the same ROAs:
 - a. Select the replicates, Click  (Define ROA) in the toolbar above the melt curve plots, then click-drag an area in one of the plots to define a melt phase and replace the ROA. Repeat for each melt phase you identify.

- b. If necessary, adjust the start and end temperatures for each ROA:
 - **To move the ROA:** Starting from within the ROA, click-drag the ROA.
 - **To adjust the start and end temperatures individually:** Click-drag the ROA start or the end line.



IMPORTANT! Make sure that the fluorescence at the start temperature is lower than the fluorescence at the end temperature.

3. Click  **Analyze** to reanalyze using the edited ROAs.

Note: After you edit ROAs, the analysis mode changes from Auto to Manual.

Review the flags and Tm values

Review the flags and Tm values, then consider editing the analysis settings and/or omitting wells before you review the replicate results.

1. In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann Tm
 - Derivative Tm
 - Unselected Wells
 - Legend
2. Review flagged wells in the Well Table:
 - a. Click the  (Flag Indicator) column header to sort the wells according to the number of flags applied to the well.
 - b. Scroll the table to the right to view the flags that are applied to the wells.



- c. For flagged wells, select the well in the Well Table, then review the melt curves for the well, compared to the other wells in the replicate group.
- d. Omit wells from analysis as necessary.


- For each replicate group, select all the wells in the replicate group, then review the Tm B (Boltzmann Tm) and the Tm D (derivative Tm) in the Well Table and in the melt curves:

Note: In the melt curves, the Boltzmann Tm is a green dashed vertical line, and the derivative Tm is black dotted vertical line.

- Are the Tm B or Tm D values significantly different from the Tm values for other wells in the replicate group? Do any replicates have melt curves that are inconsistent with the other melt curves for the replicate group? If so, you may consider omitting wells from analysis.
- Are the melt curves within the replicate group similar?

- If you omitted any wells from the analysis, click ► **Analyze**.

Review the Boltzmann fit

- In the Analysis > Well Results screen, select  **Show in Plot**, then select to show these plot components:
 - Boltzmann fit
 - Unselected Wells
 - Legend
- Scroll through each well in the Well Table, then compare the fluorescence melt curve to the Boltzmann fit curve (dark green thick curve) and review the value in the B Fit (Boltzmann fit) column of the Well Table:
 - How well does the fluorescence melt curve correspond to the Boltzmann fit curve?
 - Is the B Fit value close to 1?
 - Is the B Fit value similar among wells in the replicate group?

Note: When you define the ROA manually, you may observe a gap between the ROA start or end temperature and the start or end of the Boltzmann fit curve. The gap occurs if the defined ROA start or end temperature does not correspond exactly with a fluorescence datapoint.

Example well results

For the ViiA™ 7 System example ligand screening study, observe the following:

- Fluorescence levels are flat in the NPC and LOC wells.
- For the sample and reference replicate groups, the fluorescence levels as displayed in the fluorescence melt curve are not significantly different, so you do not need to edit the baseline.
- The sample and reference wells contain one peak in the derivative melt curve.
- Within each replicate group:
 - The fluorescence melt curves and the Boltzmann Tm values are similar.
 - The derivative melt curves and the derivative Tm values are similar and there are no outliers.
- For each well, the region of analysis meets the recommended criteria.


Review the replicate results

Using the Protein Thermal Shift™ Software, review the T_m statistics for the replicate groups and look for positive hits in the Analysis > Replicate Results screen.


This section provides guidance on how to review and interpret the replicate results. Some common troubleshooting causes are provided. For more detailed information on troubleshooting, see Chapter 8, “Troubleshooting”.

Review the T_m statistics


In the Analysis > Replicate Results screen, review the T_m statistics to evaluate the variability among replicates. The T_m statistics for each ROA are plotted along the x-axis for each replicate group.

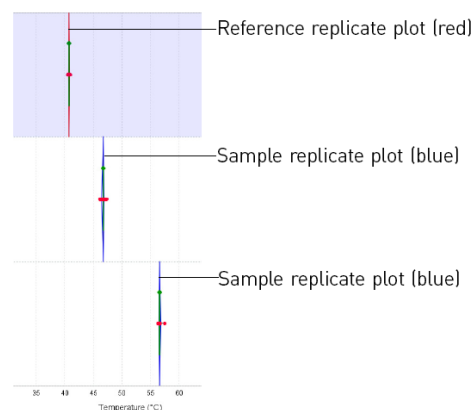
1. In the Analysis > Replicate Results screen, select  **Plot by**, then select the type of T_m statistics to review:
 - **T_m-Boltzmann**
 - **T_m-Derivative**
2. Specify the condition hierarchy to group the replicate plots and change the order of conditions in the plot:


Note: Changing the condition hierarchy does not affect the results; it only changes how the replicate plots are grouped and the order in which the conditions are displayed in the Replicate Results Plot.

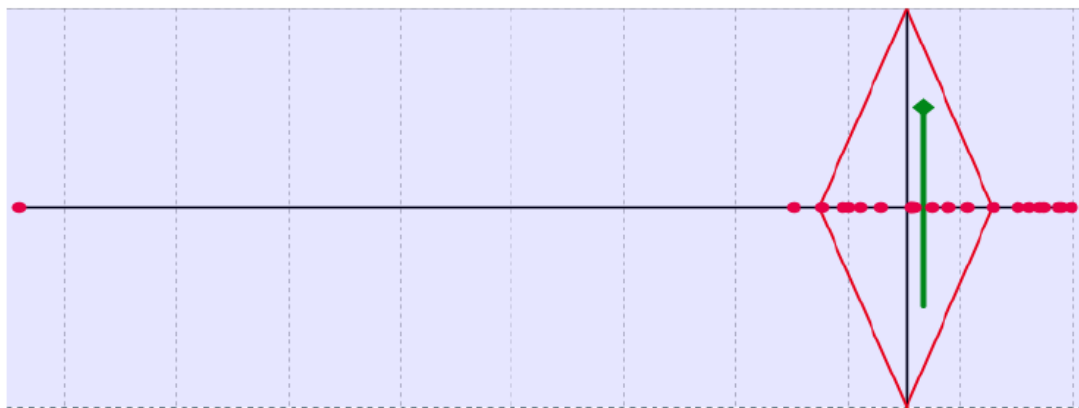
- a. Click  **Condition Hierarchy** above the top right corner of the plot.
- b. In the dialog box, select a condition, then use the Up and Down arrows to change the hierarchy of conditions.
The condition at the top-most level of the hierarchy is displayed on the far-right side of the Replicate Results Plot, and the replicate plots are grouped according to the top-most condition.

Note: For the example ligand screening study, set up the hierarchy so that **Ligand** is at the bottom in the dialog box and displayed on the far-left side of the Replicate Results Plot.

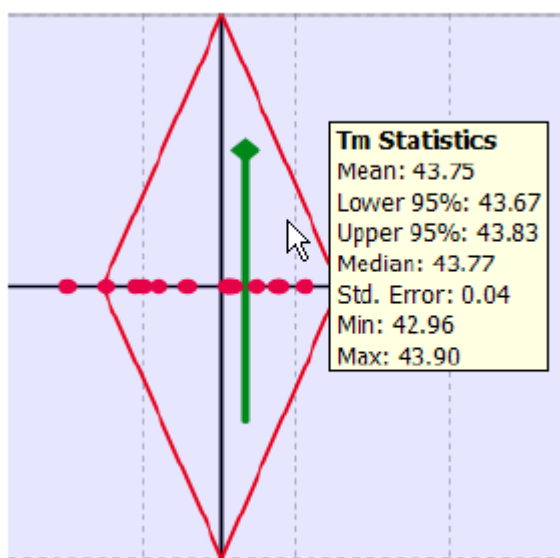
3. Scan the Replicate Results Plot to review the conditions that affect the T_m values relative to the reference replicate group.
4. Review the T_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot one or more times.



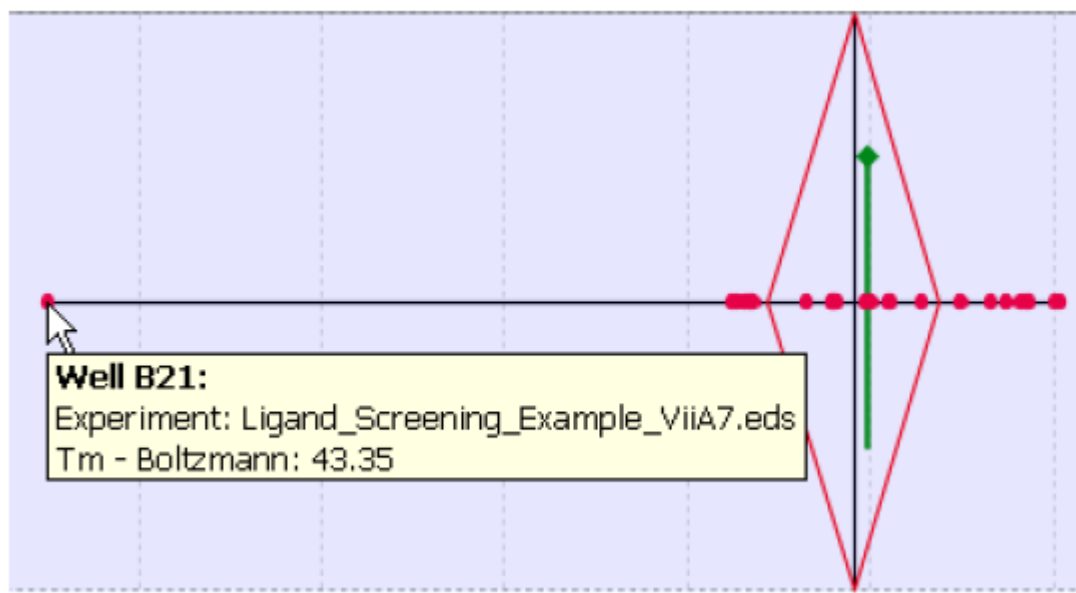
- b. Move the plot: Click , then click-drag the plot until the replicate plot of interest is in view.




- c. Place the cursor within the diamond, then wait to view a tooltip with the Tm statistics for the replicate group.



- d. To examine outliers, place the cursor over a datapoint, then wait to view a tooltip with the well information, experiment file name, and the T_m selected for the plot.



Note: Click  to restore the default zoom.


5. In the Replicate Groups table, review the T_m statistics for each replicate group:
- **Std. Error** (standard error of the mean for the T_m value): Is the value low? If the value is high, review the data for each replicate.
 - **Min** and **Max** (minimum and maximum T_m values for the replicate group): Is the range of T_m values for the replicate group within 1 degree? If the range of T_m values exceeds 1 degree, review the data for each replicate.
6. Omit outliers as necessary, then click  **Analyze**.


Review the positive hits

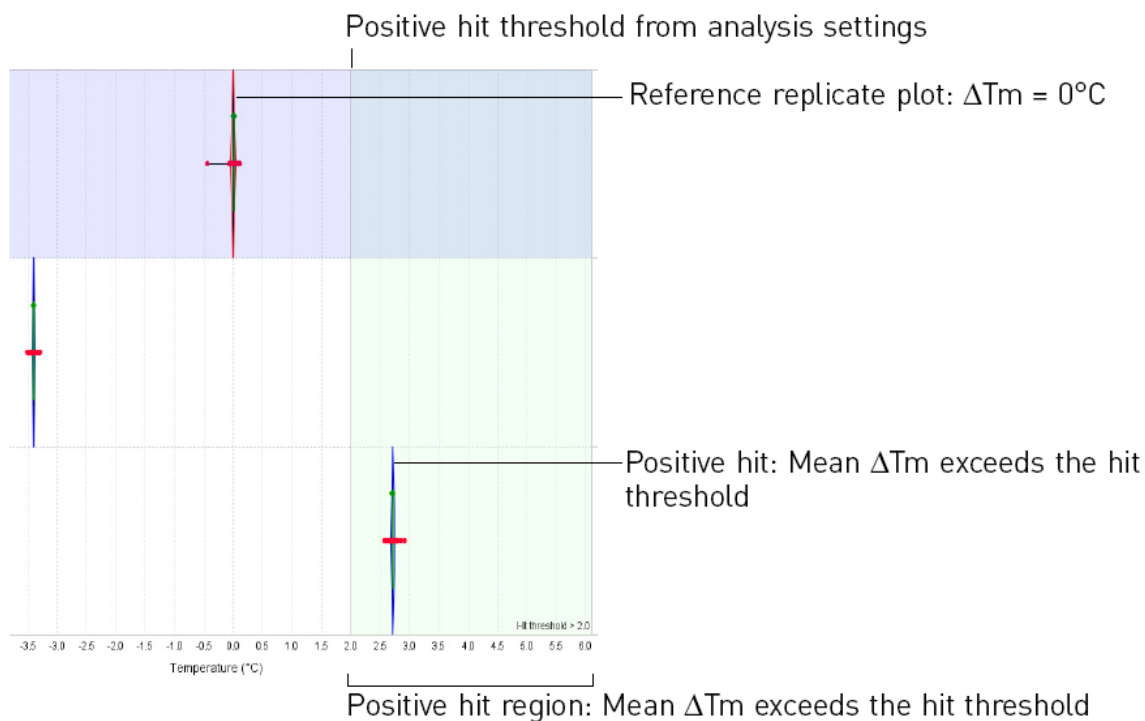
In the Analysis > Replicate Results screen, review the positive hits to identify the conditions that produce the maximum effect on thermal stability. Replicate groups with positive hits have ΔT_m values that exceed the threshold set in the analysis settings.



Note: The positive hits are determined according to the analysis settings that you specified.

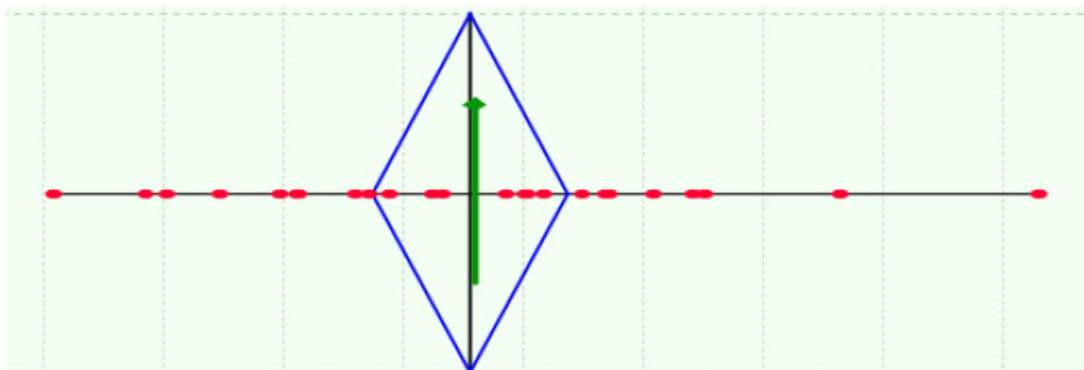
Note: You must specify the reference replicate group to calculate ΔT_m values and to determine positive hits.

1. In the Analysis > Replicate Results screen, select  **Plot by**, then select the type of ΔT_m statistics to review:
- **ΔT_m -Boltzmann**
 - **ΔT_m -Derivative**

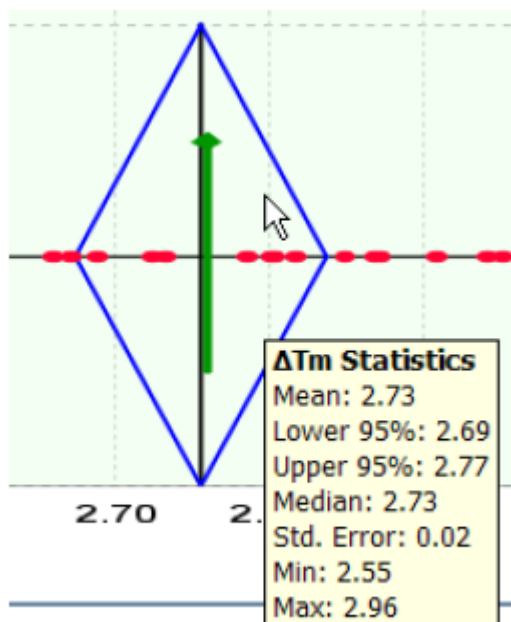
2. Select the  **Show in Plot** menu, then select to show **Positive Hits**.
3. Scan the green shaded area of the Replicate Results Plot for positive hits.



4. Review the ΔT_m statistics for each replicate group in the Replicate Results Plot:
 - a. Zoom in: Click , then click-drag an area on the plot as many times as you need to magnify the plot.
 - b. Move the plot: Click , then click-drag the plot until the replicate plot you want to review is in view.



- c. Place the cursor over the replicate plot, then wait to view a tooltip with the ΔT_m statistics for the replicate group.



Note: Click  to restore the default zoom.

5. In the Replicate Groups table, review the positive hits (✔) in the Hits B or Hits D column and review the ΔT_m B or ΔT_m D statistics in the table.



Troubleshooting

Observation	Possible cause	Recommended action
High initial background signal and/or a small transitional increase in signal	Native protein has external hydrophobic sites	The protein may not be a suitable candidate for Protein Thermal Shift™ studies. Perform protein:dye titration studies to optimize the protein concentration and protein:dye ratio.
Flat signal or decrease in signal	Protein is heat-stable and the T _m exceeds the range for Protein Thermal Shift™ studies (98°C or higher)	Use an alternate method to screen for conditions that increase thermal stability of the protein. You may consider performing Protein Thermal Shift™ studies to screen for buffers, ligands, or mutations that decrease thermal stability of the protein.
High initial background signal High fluorescence in NPC wells	Protein solution contains high levels of detergent (> 0.02%)	Perform protein:dye titration studies to optimize the protein concentration and protein:dye ratio. Repurify the protein using an ammonium sulfate precipitation method. Resolubilize the purified protein using HEPES buffer or a buffer with neutral pH, then add glycerol and DTT.
High initial background signal High fluorescence in NPC wells High fluorescence in LOC wells	Buffer component interacts with the dye	Perform protein:dye titration studies to optimize the protein concentration and protein:dye ratio. Perform a buffer screening Protein Thermal Shift™ study to identify alternative buffer conditions.
High initial background signal High fluorescence in LOC wells	The ligand interacts with the dye	Use an alternate method to screen for conditions that affect thermal stability of the protein.
High initial background signal Flat signal or decrease in signal	Protein aggregation or the protein is partially unfolded	Repeat the study with a fresh protein sample. Perform a buffer screening Protein Thermal Shift™ study to identify buffer conditions that increase thermal stability of the protein, then repeat the original study using the new buffer conditions.
Melt curves in NPC wells Melt curves in LOC wells	Protein contamination in the buffer, buffer component, ligand, or dye	Repeat the study with fresh reagents.

(continued)

Observation	Possible cause	Recommended action
Melt curves with low relative fluorescence levels	Protein concentration is too low	Perform protein:dye titration studies to optimize the protein concentration and protein:dye ratio.
Flat signal or decrease in signal	Passive reference is not set to None in the experiment file	<ol style="list-style-type: none"> Using your real-time PCR system software, open the experiment file. Set the passive reference to None. <ul style="list-style-type: none"> ViiA™ 7 Software: Define screen StepOne™ Software or 7500 Software: Plate Setup > Assign Targets and Samples screen Save and reanalyze the experiment. Reimport the experiment file into the Protein Thermal Shift™ study.
No fluorescence or very low fluorescence	Incorrect filters selected with the ViiA™ 7 Real-Time PCR System.	Repeat the protein melt reactions and instrument run, making sure that you select the correct filters when you set up the experiment file (see “Create and set up an experiment file for the ViiA™ 7 Real-Time PCR System” on page 28).

Examples of observations

High initial background signal

In this example, protein samples with high detergent levels produce a high initial background signal in the melt curves.

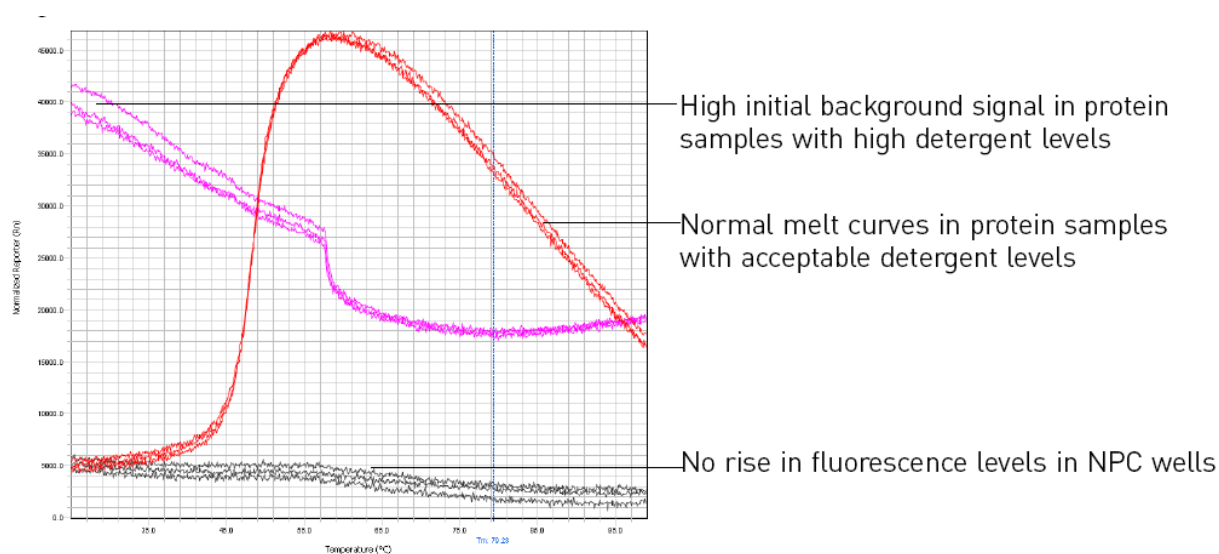


Figure 2 Melt curves in ViiA™ 7 Software

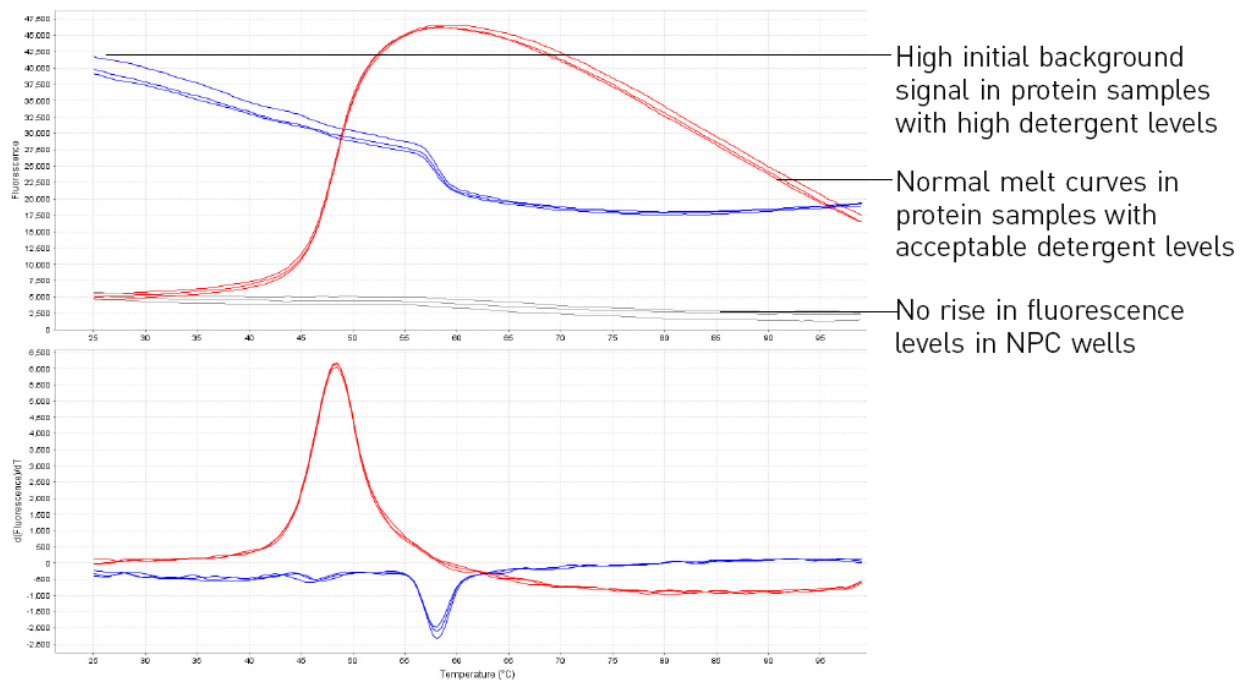
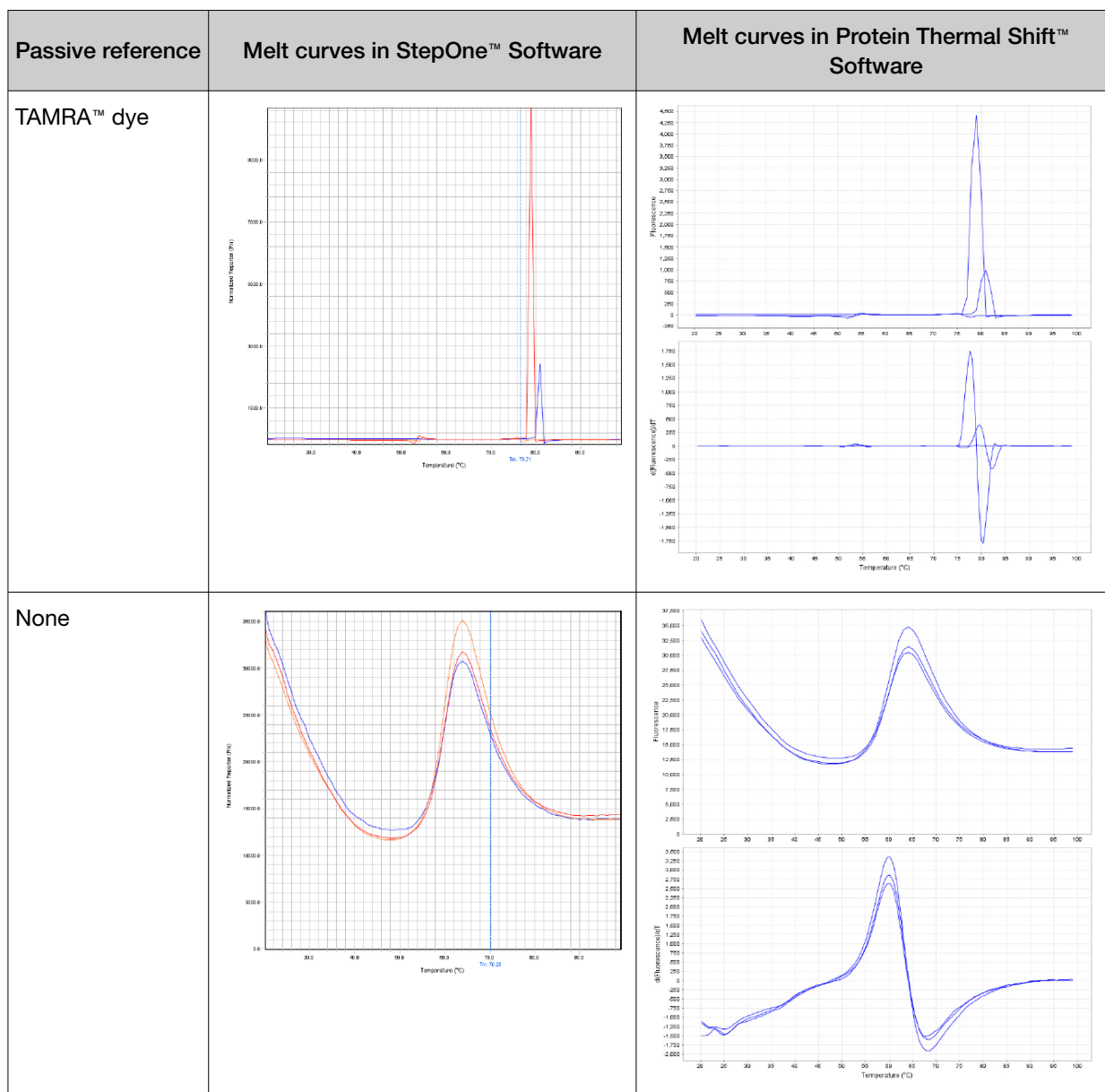


Figure 3 Melt curves in Protein Thermal Shift™ Software

Flat signal or decrease in signal

In this example, flat signals in the melt curves are observed because the passive reference is not set to None.





Supplemental information

Hardware recommendations

The Protein Thermal Shift™ Software v1.4 is fully operational when installed on a computer with the recommended hardware configuration.

Component	Recommended configuration	Minimum requirements ^[1]
Computer	<ul style="list-style-type: none">• 2.4 GHz CPU• 2GB of RAM• Disk space:<ul style="list-style-type: none">– One hard drive (no partitions): 20 GB free space– Two hard drives or two partitions: 300 MB free space on the Programs drive, 20 GB free space on the data drive, and 1 GB free space on the user files drive	<ul style="list-style-type: none">• Intel™ Pentium® 4 processor or compatible processor, 1.2 GHz• 1GB of RAM• One hard drive with 10 GB free space
Monitor	<ul style="list-style-type: none">• 1280 × 1024 pixel resolution for full screen display• 16-inch color monitor• 32-bit color	<ul style="list-style-type: none">• 1280 × 1024 pixel resolution for full screen display• 16-inch color monitor• 32-bit color
Operating system	Microsoft™ Windows™ 10 Operating System 32-bit, Service Pack 2 or later, or Microsoft™ Windows™ 10 Operating System, 32-bit	Microsoft™ Windows™ 10 Operating System 32-bit, Service Pack 2 or later

^[1] The Minimum requirements column lists the lowest specifications that permit the software installation. The minimum requirements may not provide optimal performance. The support of an installation in this environment is not guaranteed.

Boltzmann fitting method

Data from the region of analysis are fit to the Boltzmann equation to generate the T_m .

The Boltzmann equation:

$$F(T) = F(pre) + \frac{[F(post) - F(pre)]}{1 + e^{\frac{(T_m - T)}{C}}}$$

where:

- $F(T)$ is the fluorescence at a particular temperature
- $F(pre)$ is the fluorescence before the transition or melting at the start of the region of analysis (ROA)
- $F(post)$ is the fluorescence after the transition or melting at the end of the ROA
- T_m is the melting temperature
- C is the enthalpy of the reaction



Software warranty information

Computer configuration

Life Technologies™ supplies or recommends certain configurations of computer hardware, software, and peripherals for use with its instrumentation. Life Technologies™ reserves the right to decline support for or impose extra charges for supporting nonstandard computer configurations or components that have not been supplied or recommended by Life Technologies™. Life Technologies™ also reserves the right to require that computer hardware and software be restored to the standard configuration prior to providing service or technical support. For systems that have built-in computers or processing units, installing unauthorized hardware or software may void the Warranty or Service Plan.

Limited product warranty

Limited warranty

Life Technologies™ warrants that for a period of ninety (90) days from the date the warranty period begins, its Protein Thermal Shift™ Software v1.4 will perform substantially in accordance with the functions and features described in its accompanying documentation when properly installed on the instrument system for which it is designated, and that for a period of ninety (90) days from the date the warranty period begins, the tapes, diskettes, or other media bearing the software product will be free of defects in materials and workmanship under normal use. If buyer believes that it has discovered a failure of the software to satisfy the foregoing warranty, and if buyer notifies Life Technologies™ of such failure in writing during the ninety (90) day warranty period, and if Life Technologies™ is able to reliably reproduce such failure, then Life Technologies™, at its sole option, will either (i) provide any software corrections or bug-fixes of the identified failure, if and when they become commercially available, to buyer free of charge, or (ii) notify buyer that Life Technologies™ will accept a return of the software from the buyer and, upon such return and removal of the software from buyers systems, terminate the license to use the software and refund the buyers purchase price for the software. If there is a defect in the media covered by the above warranty and the media is returned to Life Technologies™ within the ninety (90) day warranty period, Life Technologies™ will replace the defective media. Life Technologies™ does not warrant that the software will meet buyers requirements or conform exactly to its documentation, or that operation of the software will be uninterrupted or error free.

Warranty period effective date

Any applicable warranty period under these sections begins on the earlier of the date of installation or ninety (90) days from the date of shipment for software installed by Life Technologies™ personnel. For all software installed by the buyer or anyone other than Life Technologies™, the applicable warranty period begins the date the software is delivered to the buyer.

Warranty claims

Warranty claims must be made within the applicable warranty period.

Warranty exceptions

The above warranties do not apply to defects resulting from misuse, neglect, or accident, including without limitation: operation outside of the environmental or use specifications, or not in conformance with the instructions for the instrument system, software, or accessories; improper or inadequate maintenance by the user; installation of software or interfacing, or use in combination with software or products, not supplied or authorized by Life Technologies™; and modification or repair of the product not authorized by Life Technologies™.

The foregoing provisions set forth Life Technologies™' sole and exclusive representations, warranties, and obligations with respect to its products, and Life Technologies™ makes no other warranty of any kind whatsoever, expressed or implied, including without limitation, warranties of merchantability and fitness for a particular purpose, whether arising from a statute or otherwise in law or from a course of dealing or usage of trade, all of which are expressly disclaimed.

Warranty limitations

The remedies provided herein are the buyer's sole and exclusive remedies. Without limiting the generality of the foregoing, in no event shall Life Technologies™ be liable, whether in contract, tort, warranty, or under any statute (including without limitation any trade practice, unfair competition, or other statute of similar import) or on any other basis, for direct, indirect, punitive, incidental, multiple, consequential, or special damages sustained by the buyer or any other person or entity, whether or not foreseeable and whether or not Life Technologies™ is advised of the possibility of such damages, including without limitation, damages arising from or related to loss of use, loss of data, failure or interruption in the operation of any equipment or software, delay in repair or replacement, or for loss of revenue or profits, loss of good will, loss of business, or other financial loss or personal injury or property damage.

No agent, employee, or representative of Life Technologies™ has any authority to modify the terms of this Limited Warranty Statement or to bind Life Technologies™ to any affirmation, representation, or warranty concerning the product that is not contained in this Limited Warranty Statement, and any such modification, affirmation, representation, or warranty made by any agent, employee, or representative of Life Technologies™ will not be binding on Life Technologies™ unless in a writing signed by an executive officer of Life Technologies™.

This warranty is limited to the buyer of the product from Life Technologies™ and is not transferable.

Some countries or jurisdictions limit the scope of or preclude limitations or exclusion of warranties, of liability, such as liability for gross negligence or willful misconduct, or of remedies or damages, as to the extent set forth above. In such countries and jurisdictions, the limitation or exclusion of warranties, liability, remedies or damages set forth above shall apply to the fullest extent permitted by law, and shall not apply to the extent prohibited by law.



Ordering information

For more information on the Protein Thermal Shift™ reagents and Protein Thermal Shift™ Software, go to [thermofisher.com](https://www.thermofisher.com).

For information to order plates and accessories, go to [thermofisher.com/plastics](https://www.thermofisher.com/plastics).

Table 1 Products for Protein Thermal Shift™ studies

Item	Cat. No.
Protein Thermal Shift™ Starter Kit	4462263
Protein Thermal Shift™ Dye Kit	4461146
Protein Thermal Shift™ Software v1.4 – Additional License	4466038
Protein Thermal Shift™ Software v1.4, 10 licenses	4466037



Glossary

analysis group

In Protein Thermal Shift™ studies, one reference replicate group and multiple sample replicate groups analyzed together to calculate ΔT_m values and to determine positive hits.

analysis mode

In the Protein Thermal Shift™ Software, the method of defining the regions of analysis (ROAs):

- **Auto:** The software defines the ROA(s):
 - **Auto: Single T_m:** The software detects one ROA and determines one T_m within the ROA.
 - **Auto: Multiple T_m:** The software detects more than one ROA and determines one T_m for each ROA.
- **Manual:** Each ROA is manually defined within the Protein Thermal Shift™ Software

The analysis mode is displayed in the Analysis Mode column of the well table in the Analysis screens.

Boltzmann fit (B fit)

In Protein Thermal Shift™ studies, a value that indicates how tightly the fluorescence data within the ROA corresponds to the Boltzmann equation.

Boltzmann T_m (T_m B)

In Protein Thermal Shift™ studies, the melting temperature (°C) calculated by fitting data in the region of analysis to the Boltzmann equation.

condition

In Protein Thermal Shift™ studies, a set of values to define a component of the protein melt reaction. For example, the default conditions in the Protein Thermal Shift™ Software are Protein, Ligand, Buffer, and Salt.

You can define the conditions and the condition values in the Setup > Conditions screen of the software.

derivative T_m (T_m D)

In Protein Thermal Shift™ studies, the melting temperature (°C) calculated for the region of analysis using the derivative of the melt curve.

experiment file

An electronic record that contains all information about a particular plate, including metadata (name, barcode, comments), plate setup (well contents, assay definitions), run method (thermal profile), run results, analysis settings, analysis results, and other plate-specific data. The analysis settings and analysis results in the experiment files are limited to the protein melt data from the real-time PCR instrument. The experiment files do not contain any settings or results from the Protein Thermal Shift™ analysis. Experiment files have the suffixes .eds (experiment document single), .edt (template), and .edm (multiple).

export

A software feature that allows you to save raw data and analysis results to *.txt or *.csv files. You can select which data to export, the file format, and the file name and location.

LOC

Ligand Only Control. Protein melt reactions that contain only ligand, buffer, water, and dye.

melt curve plot

A plot of fluorescence data collected during the melt curve stage. Melt curve plots are displayed in the real-time PCR software and the Protein Thermal Shift™ Software.

The melt curves in the real-time PCR software may not exactly match the melt curves in the Protein Thermal Shift™ Software. When the experiment files are imported into the Protein Thermal Shift™ Software, the Protein Thermal Shift™ Software reduces the noise in the fluorescence data.

In the Protein Thermal Shift™ Software, the shape of the melt curve can indicate the protein thermal stability or multiple states of protein-ligand binding. You can view the melt curve as determined by fitting the fluorescence data to the Boltzmann equation (in single T_m mode) or as a differential melt curve (in single T_m mode or in multiple T_m mode).

melt phase

In Protein Thermal Shift™ studies, a region within the melt profile that shows a nearly continuous increase in fluorescence followed by a noticeable decrease in stabilization of fluorescence levels. A melt profile may show one or more melt phases.

melting temperature (T_m)

The temperature (°C) at which 50% of the protein is folded and 50% of the protein is melted. The T_m values are displayed in the melt curve plots.

NPC

No Protein Control. Protein melt reactions that contain only buffer, water, and dye.

omit well

An action that you perform in the Protein Thermal Shift™ Software before reanalysis to omit one or more wells from calculations. Omitted wells are analyzed and can be displayed in the Replicate Results plot. The results are hidden from the results tables and are not used to calculate T_m statistics for the associated replicate group. You can add wells back to the analysis; no information is permanently discarded.

outlier

A measurement that deviates significantly from the measurement of the other replicates in the replicate group.

plate layout

An illustration of the grid of wells and assigned content in the reaction plate. The number of rows and columns in the grid corresponds to the instrument reaction block that you use.

In the Protein Thermal Shift™ Software, you can use the plate layout as a selection tool to make or view condition assignments or to omit wells from analysis. The plate layout can be saved as a *.jpg or a *.png image file.

positive hit



In Protein Thermal Shift™ studies, a sample replicate group with an average ΔT_m value that exceeds the threshold set in the analysis settings for the study. A positive hit can identify conditions (for example, buffer, ligand, or mutation) that increase or decrease protein thermal stability relative to the reference replicate group.

reference

In Protein Thermal Shift™ studies, the task assigned to one replicate group within an analysis group. The results from the reference replicate group are used to calculate ΔT_m values for sample replicate groups.

region of analysis

A range of temperatures in the melt curves from which the fluorescence data are used to calculate the T_m . The region of analysis can be detected by the software (auto analysis mode) or manually defined by the user (manual analysis mode). Up to 6 regions of analysis can be defined for a well.

replicate group

In Protein Thermal Shift™ studies, a set of identical reactions with the same set of conditions, the same task, and the same analysis group. We recommend that you include at least 4 replicates in a replicate group.

ROA

See region of analysis on page 125.

study

In the Protein Thermal Shift™ Software, a study contains one or more experiment files with one or more analysis groups. For each analysis group, the goal of the study is to identify condition sets that have the greatest ΔT_m effect on protein thermal stability relative to the reference.

task

In the Protein Thermal Shift™ Software, the type of reaction performed in the well. Available tasks in the software:

- **Sample:** Wells with protein melt conditions that you are testing
- **Reference:** Wells with protein melt conditions to use as the basis for calculating ΔT_m values and determining positive hits
- **NPC (No Protein Control):** Wells that contain no protein

T_m

See melting temperature (T_m) on page 124.

T_m B

See Boltzmann T_m (T_m B) on page 123.

T_m D

See derivative T_m (T_m D) on page 123.



Safety



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, visit thermofisher.com/support.

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the "Documentation and Support" section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if needed) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



WARNING! HAZARDOUS WASTE (from instruments). Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



WARNING! 4L Reagent and Waste Bottle Safety. Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position.

Biological hazard safety



WARNING! Potential Biohazard. Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 6th Edition, HHS Publication No. (CDC) 300859, Revised June 2020
<https://www.cdc.gov/labs/pdf/CDC-BiosafetymicrobiologicalBiomedicalLaboratories-2020-P.pdf>
- Laboratory biosafety manual, fourth edition. Geneva: World Health Organization; 2020 (Laboratory biosafety manual, fourth edition and associated monographs)
www.who.int/publications/i/item/9789240011311



Documentation and support

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

